

RESEARCH ARTICLE

# Differential immunomodulation in human monocytes versus macrophages by filarial cystatin

Gopinath Venugopal, Marion Mueller, Susanne Hartmann\*, Svenja Steinfelder

Institute of Immunology, Centre for Infection Medicine, Freie Universität Berlin, Berlin, Germany

\* [susanne.hartmann@fu-berlin.de](mailto:susanne.hartmann@fu-berlin.de)



**OPEN ACCESS**

**Citation:** Venugopal G, Mueller M, Hartmann S, Steinfelder S (2017) Differential immunomodulation in human monocytes versus macrophages by filarial cystatin. PLoS ONE 12(11): e0188138. <https://doi.org/10.1371/journal.pone.0188138>

**Editor:** Jyotika Sharma, University of North Dakota, UNITED STATES

**Received:** June 3, 2017

**Accepted:** November 1, 2017

**Published:** November 15, 2017

**Copyright:** © 2017 Venugopal et al. This is an open access article distributed under the terms of the [Creative Commons Attribution License](https://creativecommons.org/licenses/by/4.0/), which permits unrestricted use, distribution, and reproduction in any medium, provided the original author and source are credited.

**Data Availability Statement:** All relevant data are within the paper and its Supporting Information files.

**Funding:** This work was supported by the German Research Foundation (DFG - International Research Training Group 1673) and Sonnenfeld Foundation, Berlin, Germany. The funders had no role in study design, data analysis, decision to publish or preparation of the manuscript.

**Competing interests:** The authors have declared that no competing interests exist.

## Abstract

Parasitic nematodes have evolved powerful immunomodulatory molecules to enable their survival in immunocompetent hosts by subverting immune responses and minimizing pathological processes. One filarial molecule known to counteract host immune responses by inducing IL-10 and regulatory macrophages in mice is filarial cystatin. During a patent filarial infection monocytes encounter microfilariae in the blood, an event that occurs in asymptotically infected filariasis patients that are immunologically hyporeactive. The microfilarial larval stage was formerly shown to induce human regulatory monocytes and macrophages. Thus, here we aim was to determine how filarial cystatin of the human pathogenic filaria *Brugia malayi* (BmCPI-2) contributes to immune hyporesponsiveness in human monocytes and macrophages elicited by microfilaria. For this purpose, filarial cystatin was depleted from microfilarial lysate (Mf). Detecting the immunomodulatory potential of cystatin-depleted Mf revealed that IL-10, but not IL-8 and IL-6 induction in monocytes and macrophages is dependent on the presence of cystatin. In addition, the Mf-induced expression of the regulatory surface markers PD-L1 and PD-L2 in human monocytes, but not in macrophages, is dependent on cystatin. While Mf-treated monocytes result in decreased CD4<sup>+</sup> T-cell proliferation in a co-culture assay, stimulation of T-cells with human monocytes treated with cystatin-depleted Mf lead to a restoration of CD4<sup>+</sup> T-cell proliferation. Moreover, IL-10 induction by cystatin within Mf was dependent on p38 and ERK in macrophages, but independent of the ERK pathway in monocytes. These findings indicate that filarial nematodes differentially trigger and exploit various signaling pathways to induce immunomodulation in different myeloid cell subsets.

## Introduction

Helminths play a central role in causing infections among communities living in tropical and subtropical regions, and about one third of the world population is infected with one or more parasitic helminths[1]. The two major classes of helminths are nematodes and platyhelminths, both of which contain causative agents of a variety of acute debilitating diseases and chronic conditions[2]. In general, helminth infections are well characterized by strong Th2 immune responses, and are able to down-regulate host immune responses[1,3,4].

Lymphatic filariasis (LF) is one of seventeen Neglected Tropical Diseases (NTDs) classified by the World Health Organization (WHO)[5], and is the third leading parasitic cause of morbidity and disability worldwide. LF is normally associated with varied clinical outcomes and according to WHO Global Health Estimates (GHE), in 2015 alone LF contributed close to 2 million Disease Adjusted Life Years (DALYs)[6]. The causative agents of human LF are three species of parasitic nematodes, namely *Wuchereria bancrofti*, *Brugia malayi*, and *Brugia timori*. Over 120 million people around the world are infected with either of these parasites and about 40 million people suffer from disfigurement and incapacitation for work as a result of LF-induced symptoms like lymphedema, elephantiasis or hydrocele[7]. The lifecycle of these filarial worm species comprises of both mosquito (*Aedes*, *Anopheles*, *Culex* or *Mansonia* spp.) and human stages. Infection with the parasite is initiated when an infected mosquito deposits the infective larvae stage on the skin of a human host during a blood meal. The parasite larvae then reach the afferent lymphatic vessels, where they are able to undergo two moulting stages and eventually mature to adult worms. After mating, females start producing microfilariae, which are then able to circulate in the peripheral blood of the human host. During a subsequent blood meal the mosquito vector is therefore able to pick up circulating blood microfilariae, thus initiating the development of the infective L3 stage larvae and completing the transmission cycle of the parasite.

The host-parasite interaction in LF often results in a chronic infection associated with a wide clinical spectrum, characterized by functional dysregulation of both the innate and the adaptive immune responses[8]. Three of the key clinical manifestations of this disease include asymptomatic infection, despite the presence of the larval blood stages (microfilariae) caused by hyporesponsiveness towards the parasite, chronic pathology with hydrocele, or elephantiasis, with or without active infection displaying a Th1/Th17-biased immune response, and endemic normality, where the host is putatively immune against the parasite.

The longevity of helminth parasites in the mammalian host is primarily determined by their capacity to modulate the immune system. Immunomodulatory helminth-derived products such as protease inhibitors[9–14] and homologues of mammalian cytokines and chemokines[15–18] induce a modified Th2-type immune response, which creates a beneficial microenvironment for the parasite[19,20]. Cystatins, a family of cysteine protease inhibitors, are known to regulate cysteine protease functions such as protein catabolism, antigen processing, inflammation, dystrophy and metastasis in viruses, bacteria, protozoa, fungi, plants and mammals [21–24]. Helminth cystatins are highly expressed protease inhibitors and they have been shown to possess strong immunomodulatory properties. They have a dual function, due to their ability to inhibit both parasite and host cysteine proteases. Cystatins are reversible competitive inhibitors of cysteine proteases and control protein processing and turnover, thereby inhibiting antigen processing[10,11], leading to diminished T cell priming. Secondly, they induce the immunosuppressive cytokine IL-10[9] and nitric oxide-producing regulatory macrophages[25], which can result in the inhibition of T cell proliferation, as shown in previous studies on cystatins from the filarial rodent parasites *Acanthocheilonema viteae* [9] and *Litomosoides sigmodontis* [14], and the human filarial nematode *Onchocerca volvulus* [12].

We could show in previous studies that monocytes and macrophages are target cells of filarial cystatins *in vitro* and *in vivo*[12,13]. Moreover, monocytes from asymptomatic individuals harbouring microfilariae from *Brugia malayi* upregulate programmed-death ligand 1 (PD-L1), IL-10 and IL-8, which is recapitulated in experiments using human monocytes and macrophages from non-endemic healthy donors stimulated with microfilarial lysate (Mf) *in vitro*. Importantly, macrophages differentiated in the presence of Mf lysate respond poorly to LPS, and Mf-stimulated monocytes are able to suppress CD4<sup>+</sup> T cell functions[26]. This is in line with the finding that patent lymphatic filariasis patients have impaired monocyte and CD4<sup>+</sup> T cell function, reflected by their inability to produce inflammatory cytokines in response to

activating stimuli[27,28], while elevated levels of IL-10 are observed in asymptomatic patients [26,29]. The anti-inflammatory cytokine IL-10 plays a major role in regulating inflammatory diseases such as allergies and autoimmune disorders[30], while IL-8 is an important chemokine with a key role in the pathogenesis of filarial diseases[31] and it displays elevated levels in patients with and without circulating microfilariae[32].

There are three known cystatins from the human filarial nematode *Brugia malayi* (BmCPI-1, -2 and -3). Among these three types, BmCPI-2 is expressed both in the vector and in the mammalian life stages while BmCPI-1 and -3 are exclusively expressed in the mosquito stages of the parasite [10,33]. Most importantly, BmCPI-2 is released in greater amounts by *Brugia malayi* microfilaria compared to adults stages[34]. This prompted us to investigate the immunomodulatory potential of BmCPI-2 on human monocytes and macrophages. Furthermore, we tried to determine the contribution of *Brugia malayi* cystatin in Mf-induced immunomodulation of human monocytes and macrophages, and to further demonstrate the key host signalling events used by BmCPI-2 and Mf to modulate these host cell types. In the present study, we demonstrate that the expression of IL-10 by Mf-stimulated monocytes and macrophages is dependent on cystatins. Moreover, we show that IL-10 expression in monocytes, but not in macrophages, is independent of ERK signalling, suggesting different pathways of IL-10 induction in these cell populations.

## Materials and methods

### Human blood samples

Human blood cells were derived from the German Red Cross and all experiments were approved by the ethics committee of the Charite University, Berlin, Germany.

### Filarial lysate preparation

Live *Brugia malayi* microfilariae were kindly donated by the NIAID/NIH Filariasis Research Reagent Resource Center in Athens, Georgia, USA. The worms were initially washed twice in RPMI medium containing 1% glucose, 200 U/ml penicillin and 200 µg/ml streptomycin, and incubated with the same medium at 37°C and 5% CO<sub>2</sub> for 24h, followed by washing in phosphate buffered saline (PBS). To prepare the filarial lysate, microfilariae worms were homogenised directly in a glass homogeniser and ultrasonicated on ice at an intensity of 10% for 3 min. The homogenate was centrifuged at 12,000 rpm at 4°C for 10 min and sterile filtered through a 0.22 µm filter. Protein concentration from the filtered homogenate was determined by using the Pierce™ BCA protein assay kit (Thermo Scientific, Rockford, IL, USA), as per the manufacturer's guidelines.

### Immunoprecipitation of *Brugia malayi* cystatin

An immunoprecipitation technique was performed to deplete *Brugia malayi* cystatin from protein mixtures of microfilarial lysate (Mf) by using rabbit polyclonal antibodies raised against BmCPI-2, recombinantly expressed in *E. coli*. The assay was performed using an immunoprecipitation kit from Novex (Thermo Fisher Scientific, Waltham, MA, USA), as per the manufacturer's guidelines. Briefly, 50 µl of protein G-coupled Dynabeads® were added to the buffer containing anti-BmCPI-2 antibodies (1:1000). After 10 min incubation, the supernatant was removed and beads were washed twice. Then, 100–150 µl of Mf was added to the beads. After 10–15 minutes incubation, the beads were removed, cystatin-depleted Mf was collected and the protein concentration was determined using the Pierce™ BCA protein assay kit (Thermo Scientific, Rockford, IL, USA), as per the manufacturer's guidelines.

## Cloning of the BmCPI-2 sequence into the LEXSY expression system and purification of BmCPI-2

The nucleotide sequence of BmCPI-2 was amplified from *Brugia malayi* microfilaria cDNA using specific primers (forward: 5' -gTCgACgCTTTgATTCATCgACgAg-3', reverse: 5' -ggTACCTACTgACgAgAgTACCTTTg-3') including restriction sites for cloning the sequence into the pLEXSY-sat2 plasmid of the LEXSYcon2 Expression Kit (Jena Bioscience GmbH, Jena, Germany). The amplified sequence did not include the coding region for the specific signal peptide of BmCPI-2. Cloning procedures followed the manufacturer's instructions and resulted in a monoclonal LEXSY cell strain expressing and secreting BmCPI-2 with a C-terminal hexa-histidine tag into the culture medium. Purification of the protein was performed via affinity chromatography using HisTrap™ excel columns and the ÄKTA™ pure chromatography system (GE Healthcare Bio-Science AB, Uppsala, Sweden) with a non-denaturing protocol, with imidazole as a competitive eluent. The purified protein was dialysed against PBS and sterile filtered. The protein concentration was determined using the Pierce™ BCA Protein Assay Kit (Pierce Biotechnology, Rockford, USA).

## Immunoblot analysis

SDS-PAGE was performed in 12% separating gel with 6% stacking gel. 15 µg of protein from each sample was applied per lane. Proteins were transferred to a nitrocellulose membrane for immunodetection. After blocking with 5% non-fat dry milk in TBS-Tween, membranes were incubated with rabbit anti-BmCPI-2 (1:50,000), followed by incubation with horseradish peroxidase conjugated anti-rabbit IgG (1:15,000) (Cell signalling, Danvers, MA, USA). Signals were detected by a chemiluminescence reaction using the ECL kit (Amersham, GE Healthcare Europe, GmBH). All bands were visualized by an enhanced chemiluminescence system (PEQLAB, Erlangen, Germany).

## Protein detection by silver staining

Silver staining was performed to detect proteins separated in SDS-PAGE by using Thermo Scientific Pierce Silver Stain kit (Thermo scientific, Rockford, IL, USA) using the manufacturer's protocol. Briefly, 15 µg of protein from each sample was applied per lane in SDS-PAGE. After running the gel for 1 h, the gel was washed twice with ultrapure water and then fixed twice by adding 30% ethanol and 10% acetic acid for 15 min each. After 30 min of fixation, the gel was washed twice, first with 10% ethanol then with ultrapure water. After the washing step, the gel was sensitized quickly for 1 min using silver stain sensitizer then washed twice with ultrapure water. The gel was then incubated for 30 min with a silver stain enhancer and washed briefly with ultrapure water. The gel was developed using silver stain developer solution and the development procedure was stopped immediately after the desired band intensity was achieved using 5% acetic acid. After the addition of stop solution, the gel was washed briefly with ultrapure water and incubated for 10 min at room temperature.

## Isolation of CD14<sup>+</sup> monocytes and differentiation protocol for monocyte-derived macrophages from human PBMCs

Peripheral blood mononuclear cells (PBMCs) were isolated from buffy coats by density gradient centrifugation using Pancoll human (PAN Biotech, Aidenbach, Bayern, Germany). To positively select for CD14<sup>+</sup> monocytes, 200 µl of anti-CD14 beads (Miltenyi Biotec, Bergisch-Gladbach, Germany) were added to 5-10x10<sup>8</sup> PBMCs for 20 min and cells were separated by an autoMACS classic (Miltenyi Biotec, Bergisch-Gladbach, Germany) using the program

'Possel'. Macrophages were generated *in vitro* by culturing CD14<sup>+</sup> monocytes in complete RPMI containing 10 ng/ml human M-CSF in 6-well cell culture plates at a cell concentration of  $0.33 \times 10^6$ /ml and a density of  $0.1 \times 10^6$ /cm<sup>2</sup> for 6 days, at 37°C and 5% CO<sub>2</sub>.

### Cell culture and stimulation of monocytes and macrophages

The CD14<sup>+</sup> monocytes and monocyte-derived macrophages were stimulated with parasite material or recombinant BmCPI-2 and kept for 24 h at 37°C and 5% CO<sub>2</sub>. The supernatant was collected after 24 h and stored at -20°C for further analysis. In experiments including inhibitors, the inhibitors were added to the cell culture 60 min before stimulation. The following inhibitors were used in our study: p38 inhibitor (SB203580), and MEK1/2 inhibitor (U0126) (both from Calbiochem, Merck, Darmstadt, Germany).

### Flow cytometry

Monocytes and macrophages were analysed for surface expression of PD-L1 and PD-L2. Cells were treated with FcR blocking reagent (Miltenyi Biotec, Bergisch-Gladbach, Germany), stained with fixable viability dye eFluor 780 (eBiosciences, San Diego, USA) and additionally stained for the following markers: anti-274-PE (clone MIH1), anti-273-PE (clone MIH18). Fixed cells were acquired using the FACS Canto II (BD, Franklin Lakes, USA) and analysed using FlowJo, version 10.2 (Tree Star, Ashland USA).

### Co-culture of CD4<sup>+</sup> T cells and monocytes

To perform the co-culture experiments for a functional assay of T cell proliferation, CD4<sup>+</sup> T cells were isolated from PBMCs using the CD4<sup>+</sup> T cell microbeads (Miltenyi Biotec, Bergisch-Gladbach, Germany) according to the manufacturer's instructions and sorted on an auto-MACS classic (Miltenyi Biotec, Bergisch-Gladbach, Germany) using the program 'Possel'. The autologous CD4<sup>+</sup> T cells were stained for CFSE and rested in complete RPMI for 24 h at 37°C and 5% CO<sub>2</sub>.

Monocytes were left unstimulated or stimulated with 20 µg/ml of Mf or cystatin-depleted Mf lysate for 24 h in a 5% CO<sub>2</sub> incubator at 37°C in 24 well plates.  $5 \times 10^5$  CFSE-labelled CD4<sup>+</sup> T cells were co-cultured with  $1 \times 10^5$  stimulated monocytes in 96 well flat bottom plates in the presence of 2 µg/ml soluble anti-CD3 (OKT3, eBioscience, San Diego, USA). After 5 days, the cells were washed and stained with fixable viability dye eFluor 780 (eBioscience, San Diego, USA) and anti-CD4-PerCP (clone RPA-T4, BioLegend, San Diego, USA). Fixed cells were acquired using the FACS Canto II (BD, Franklin Lakes, USA) and analysed using FlowJo, version 10.2 (Tree Star, Ashland USA).

### Cytokine detection by ELISA

IL-10, IL-8, IL-6 and TNFα proteins were measured using commercial enzyme-linked immunosorbent assay (ELISA) kits (eBioscience, San Diego, CA, USA). IL-12p40 was also measured using an ELISA kit (BioLegend, San Diego, CA, USA). All samples were measured in duplicates. Absorbance was read at 450nm with background wavelength subtracted at 570nm using a Synergy HT plate reader (BioTek, Winooski, VT, USA).

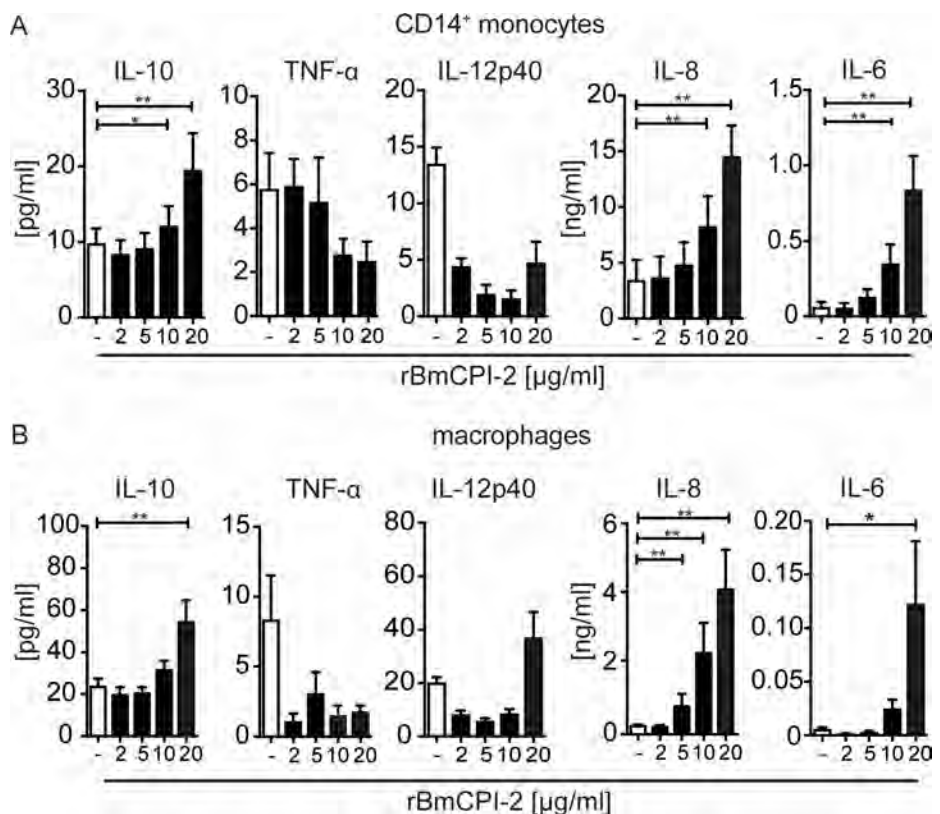
### Statistical analysis

All statistical analyses were performed using GraphPad Prism version 7.0 (GraphPad Software, Inc., CA, USA). Comparisons between groups were performed using the Wilcoxon signed rank test.

## Results

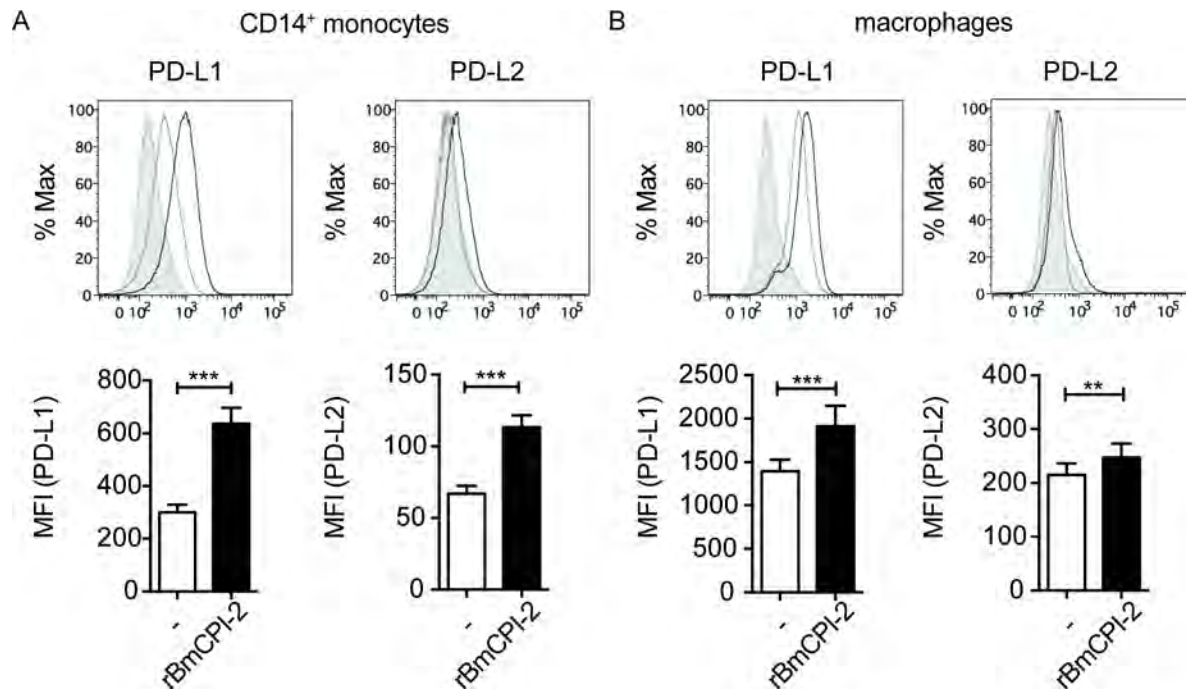
### BmCPI-2 from *Brugia malayi* induces IL-10, IL-8, IL-6 and PD-L1/L2

We have previously shown that *Brugia malayi* microfilarial lysate (Mf) induces a regulatory subset of human monocytes expressing IL-10 and PD-L1 leading to inhibition of CD4<sup>+</sup> T-cell proliferation and cytokine secretion [26]. In the present study, we hypothesized that the observed immunomodulation seen might be due to the presence of *Brugia malayi* cystatin, as the immunomodulatory potential of cystatin has been demonstrated in murine macrophages [9,14]. First, to determine the potential of BmCPI-2 to induce a regulatory phenotype in human monocytes and macrophages, we tested if BmCPI-2 is able to induce inflammatory and anti-inflammatory cytokines and programmed death-ligands in human monocytes and macrophages of healthy blood donors in a similar fashion as Mf [26]. To this end, we used recombinant BmCPI-2 (rBmCPI-2) that was expressed in a eukaryotic protein expression system. rBmCPI-2 stimulated monocytes to secrete IL-10, IL-8 and IL-6 in a dose-dependent manner, while IL-12p40 and TNF $\alpha$  were not induced (Fig 1A). Likewise, rBmCPI-2 induced IL-10, IL-8 and IL-6 in a similar fashion in human macrophages (Fig 1B). The observed increase in IL-12p40 production in human macrophages was not significant between unstimulated cells and cells stimulated with 20  $\mu$ g/ml of rBmCPI-2 (Fig 1B). Moreover, rBmCPI-2-stimulated monocytes and macrophages showed significant up-regulation of PD-L1 and PD-L2 (Fig 2).



**Fig 1. *Brugia malayi* cystatin-2 (rBmCPI-2) induces IL-10, IL-8 and IL-6 in human monocytes and macrophages.** Human monocytes (A) and macrophages (B) were stimulated with different concentrations [ $\mu$ g/ml] of recombinant BmCPI-2 for 24h. Cytokines in the supernatant were detected by ELISA. Data are from 12 donors and shown as mean  $\pm$  SEM. Statistical analysis was done using the Wilcoxon matched-pairs signed rank test in reference to the unstimulated control. \* $p$ <0.05, \*\* $p$ <0.01.

<https://doi.org/10.1371/journal.pone.0188138.g001>



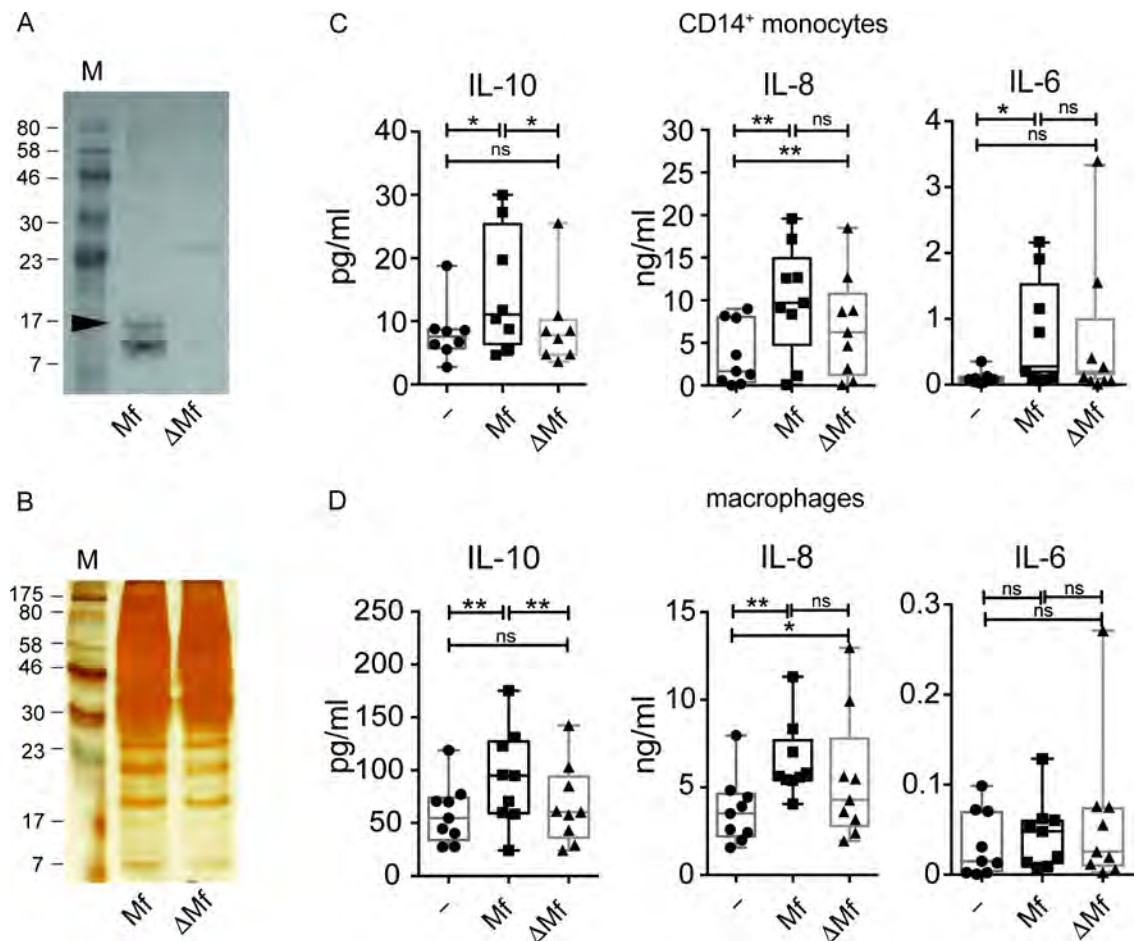
**Fig 2. rBmCPI-2 induces expression of PD-L1 and PD-L2 on human monocytes and macrophages.** Monocytes (A) or macrophages (B) were left unstimulated or stimulated for 24 h with 20  $\mu$ g/ml of recombinant BmCPI-2 (n = 12). Surface expressions of PD-L1 and PD-L2 were measured by flow cytometry. Histograms in the upper panel show representative plots of unstained (grey tinted), unstimulated (grey line) and stimulated cells (black line). Bar graphs in the lower panel show the mean fluorescence intensity (MFI). All data are represented as mean  $\pm$  SEM. P values were calculated using Wilcoxon signed-rank test. \*\*p<0.01, \*\*\*p<0.001.

<https://doi.org/10.1371/journal.pone.0188138.g002>

### Induction of IL-10 and PD-L1 and PD-L2 in human monocytes by *Brugia malayi* microfilarial lysate is dependent on cystatins

To determine if IL-10 production by microfilarial lysate (Mf) in human monocytes and macrophages is dependent on the presence of filarial cystatin, we depleted all three forms of cystatins from Mf (termed here  $\Delta$ Mf) by immunoprecipitation using polyclonal antibodies raised against recombinant BmCPI-2. BmCPI-2 was indicated by a black arrow in Fig 3A. Successful depletion of *Brugia malayi* cystatins from Mf was obtained without altering the overall protein composition of Mf (Fig 3A and 3B). Monocytes and macrophages stimulated with  $\Delta$ Mf did no longer produce significant amounts of IL-10, similar to unstimulated controls, in contrast to non-cystatin-depleted-Mf-stimulated monocytes (Fig 3C) and macrophages (Fig 3D), which induced significant amounts of IL-10. These data suggest that IL-10 induction by Mf is strictly cystatin-dependent in monocytes and macrophages.

On the other hand, Mf-induced IL-8 was not strictly dependent on cystatins, since IL-8 levels were lower but not significantly different in  $\Delta$ Mf- versus Mf-stimulated monocytes and macrophages. Similarly, Mf-induced IL-6 in monocytes was not dependent on cystatins, whereas in macrophages Mf did not induce IL-6 (Fig 3C and 3D). Notably, our results show that Mf-induced expression of PD-L1 and PD-L2 in monocytes is partially dependent on cystatins (Fig 4A). On the other hand, expression of PD-L1 in human macrophages was not dependent on cystatins (Fig 4B). Next, we employed a functional assay for CD4<sup>+</sup> T-cell proliferation using autologous monocytes treated with Mf or  $\Delta$ Mf and co-cultured both cell types with anti-CD3. We observed a significant reduction of proliferation when T-cells were co-



**Fig 3. *Brugia malayi* microfilaria lysate induced IL-10 is dependent on the presence of filarial cystatins.** Western Blot (A) and Silver gel (B) showing *Brugia malayi* microfilaria lysate (Mf) and microfilaria lysate depleted of filarial cystatins ( $\Delta$ Mf). The black arrow indicates BmCPI-2. C) Human monocytes, and D) monocyte-derived macrophages were stimulated with 20  $\mu$ g/ml of Mf or  $\Delta$ Mf for 24 h. Cytokines in the supernatant were detected by ELISA. Data are from 8 $\pm$ 9 donors and shown as mean  $\pm$ SEM. Statistical analysis was done using the Wilcoxon matched-pairs signed rank test. \* $p$ <0.05, \*\* $p$ <0.01.

<https://doi.org/10.1371/journal.pone.0188138.g003>

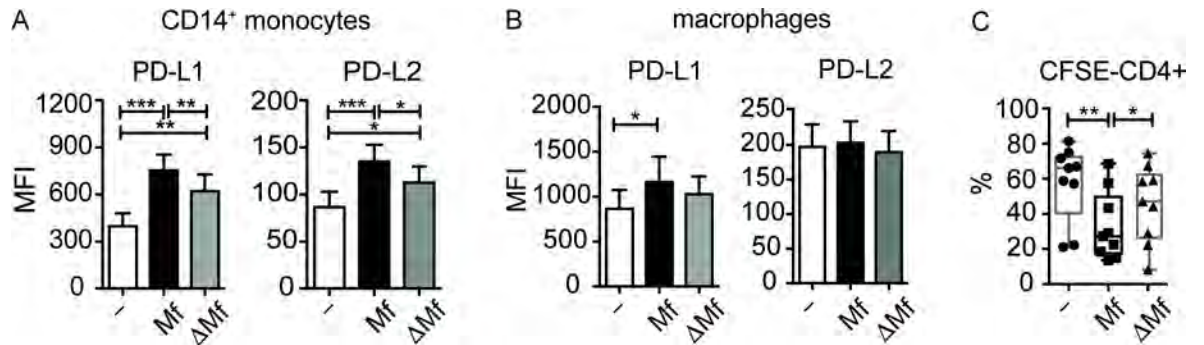
cultured with Mf-treated monocytes compared to unstimulated monocytes as described before [26]. Strikingly, when T-cells were co-cultured with  $\Delta$ Mf-stimulated monocytes proliferation was not significantly altered with respect to co-culture with unstimulated monocytes. Moreover,  $\Delta$ Mf-stimulated monocytes significantly restored T-cell proliferation compared to Mf-stimulated monocytes (Fig 4C).

Hence, IL-10 induction by Mf, but not IL-8 or IL-6, appears dependent on cystatins in both myeloid cell types. Furthermore, PD-L1 and PD-L2 expression of monocytes but not macrophages is dependent on cystatins. Most importantly, cystatins mediate the inhibitory potential of monocytes, which leads to reduced CD4<sup>+</sup> T-cell proliferation.

### Differential induction of IL-10 by rBmCPI-2-stimulated myeloid cell population

Previous studies indicate that the increased production of IL-10 in murine macrophages by cystatin from the rodent filarial nematode *Acanthocheilonema viteae* was critically dependent on two mitogen-activated protein kinases, ERK and p38[35]. Nevertheless, the immune



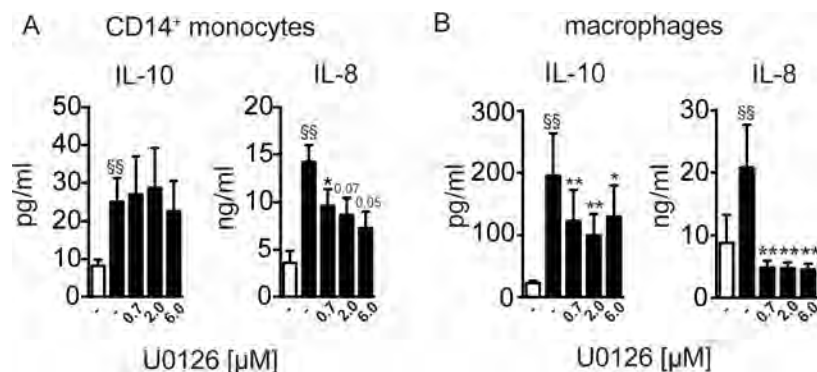


**Fig 4. Mf-induced PD-L1 and PD-L2 on monocytes is dependent on the presence of filarial cystatins.** A) Human monocytes, and B) monocyte-derived macrophages were stimulated with 20 μg/ml of Mf or ΔMf for 24 h. Surface expression of PD-L1 and PD-L2 was measured by flow cytometry. Bar graphs show the mean fluorescence intensity (MFI). Data are from 16 (monocytes) and 9 (macrophages) donors, respectively, and shown as mean ±SEM. C) Flow cytometric analysis of CD4<sup>+</sup> T cells. 5x10<sup>5</sup> CFSE-labelled CD4<sup>+</sup> T cells were co-cultured for 5 days with 1x10<sup>5</sup> monocytes left either unstimulated or stimulated for 24 h with 20 μg/ml Mf or 20 μg/ml ΔMf. Graph shows the percentage of CD4<sup>+</sup> T cells that divided. Data are from 9 donors and shown as mean ±SEM. Statistical analysis was done using the Wilcoxon matched-pairs signed rank test. \*p<0.05, \*\*p<0.01, \*\*\*p<0.001.

<https://doi.org/10.1371/journal.pone.0188138.g004>

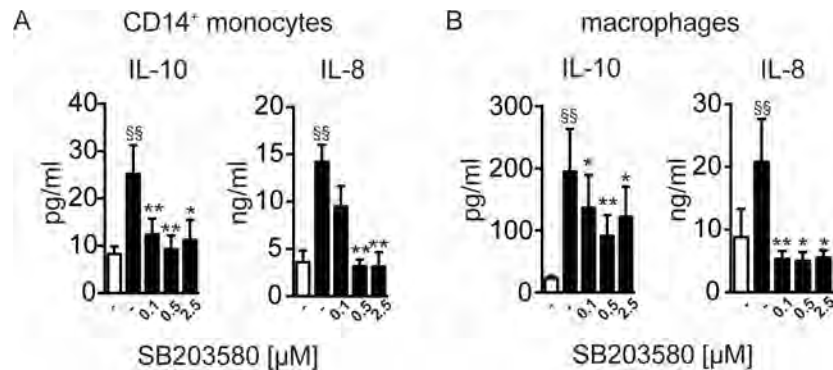
signalling events exploited by human filarial cystatin to induce IL-10 in human monocytes and macrophages is still elusive. This directed us to investigate the signalling events activated or altered by rBmCPI-2 and Mf, and leading to IL-10 and IL-8 production in human monocytes and macrophages. To this end, we used various concentrations of inhibitor molecules known to be involved in the regulation of IL-10 production[35].

Firstly, we determined the role of the MEK-dependent signalling pathway to induce IL-10 and IL-8 in rBmCPI-2-stimulated human monocytes and macrophages. Monocytes (Fig 5A) and macrophages (Fig 5B) were incubated with rBmCPI-2 for 24 h in the presence of various concentrations U0126, which is a specific inhibitor of MAP kinase kinases, MEK1 and MEK2, and thus ERK activation. Notably, rBmCPI-2-induced-IL-10 production was independent of ERK signalling, as MEK inhibition did not alter the production of IL-10 in human monocytes (Fig 5A), a finding also seen with Mf-stimulated monocytes (S1 Fig). In contrast, rBmCPI-2-induced IL-10 production in macrophages was dependent on ERK signalling, as we have observed a significant reduction of IL-10 levels in MEK inhibition (Fig 5B). Moreover, we



**Fig 5. Induction of IL-10, but not IL-8 by rBmCPI-2 in human monocytes is independent of ERK.** Human monocytes (A) and macrophages (B) were stimulated with 20 μg/ml of rBmCPI-2 for 24 h with various concentrations of MEK1/2 inhibitor. Cytokines in the supernatants were detected by ELISA. Data are from 7±9 donors and shown as mean ±SEM. Statistical analysis was done using the Wilcoxon matched-pairs signed rank test. Values statistically different from the unstimulated control are depicted as § p<0.05 and §§ p<0.01. Values statistically different from rBmCPI-2-stimulated monocytes are depicted as \* p<0.05 and \*\*p<0.01.

<https://doi.org/10.1371/journal.pone.0188138.g005>



**Fig 6. Induction of IL-10 and IL-8 by rBmCPI-2 in human monocytes and macrophages is dependent on p38.** Human monocytes (A) and macrophages (B) were stimulated with 20 μg/ml of rBmCPI-2 for 24 h with various concentrations of p38 inhibitor. Cytokines in the supernatants were detected by ELISA. Data are from 7±9 donors and shown as mean ±SEM. Statistical analysis was done using the Wilcoxon matched-pairs signed rank test. Values statistically different from the unstimulated control are depicted as § p<0.05 and §§ p<0.01. Values statistically different from BmCPI-2-stimulated monocytes are depicted as \* p<0.05 and \*\*p<0.01.

<https://doi.org/10.1371/journal.pone.0188138.g006>

observed a significant dose-dependent reduction of IL-8 levels in rBmCPI-2 (Fig 5A) stimulated monocytes and macrophages in the presence of the ERK inhibitor MEK1/2. This finding suggests that IL-10 induction by rBmCPI-2 and Mf is differentially regulated in human monocytes and macrophages.

### Induction of IL-8 and IL-10 by rBmCPI-2-stimulated myeloid cell population is dependent on p38

We next investigated the possible involvement of the p38-signalling pathway triggered by rBmCPI-2 in inducing IL-10 and IL-8 production in both human monocytes and macrophages. Inhibition of signalling pathways dependent on p38 by addition of the inhibitor SB203580 resulted in a significant reduction of IL-10 and IL-8 in both rBmCPI-2 and Mf-stimulated monocytes and macrophages (Fig 6A and 6B, S2 Fig).

## Discussion

The contribution of helminth-derived products towards the modulation of host immune responses has been addressed by several studies, and cystatin is one such secreted protein of filarial nematodes with strong immunomodulatory capacities [19,36]. With the present study we aimed to show the ability of *Brugia malayi* filarial cystatin to induce human monocytes and macrophages with a regulatory phenotype and function.

First, we demonstrated that the nematode immunomodulatory protein BmCPI-2 can significantly induce IL-10, IL-8 and IL-6, but not TNFα and IL-12p40 in human monocytes (Fig 1A) and macrophages (Fig 1B). Additionally, BmCPI-2 induced the expression of PD-L1 and PD-L2 on the surface of human monocytes and macrophages, suggesting a regulatory monocyte and macrophage phenotype *in vitro* (Fig 2). This is in line with our published data on human monocytes and macrophages stimulated with whole lysate of *Brugia malayi* microfilaria (Mf), where monocytes and macrophages develop a characteristic regulatory phenotype with high expression of IL-10 and PD-L1/L2, and inhibit CD4<sup>+</sup> T cell function [26]. Therefore, we then tested whether the observed immunomodulation seen in human monocytes and macrophages by Mf is due to the presence of cystatin. To that end we analysed the modulatory capacity of Mf that was depleted of cystatin (ΔMf). Interestingly, using polyclonal serum raised

against recombinantly expressed BmCPI-2, the protocol resulted in the depletion of all cystatins formerly detectable in Mf. Our results show that Mf-induced IL-10, but not IL-8 and IL-6 production in monocytes and macrophages is due to the presence of cystatin (Fig 3C and 3D). This is in agreement with previous studies on the ability of filarial cystatins to induce IL-10 production in murine macrophages[9,37] and human monocytes[12]. Additionally, expression of PD-L1 and PD-L2 in monocytes but not in macrophages is dependent on the presence of cystatin in Mf. Finally, both, the reduced expression of IL-10 and PD-L1/2 in monocytes, which is observed in Mf depleted of cystatins, might contribute to the observed restoration of CD4<sup>+</sup> T-cell proliferation (Fig 4).

Recently, many studies have been published addressing the therapeutic potential of cystatin in inflammatory diseases due to its anti-inflammatory activity. The modulation of innate immune cells and the profound suppression of the host immune system due to a characteristic anti-inflammatory milieu with increased levels of IL-10 and transforming growth factor (TGF)- $\beta$  in chronic helminth infections can protect against allergic diseases[38–40]. The importance of IL-10 and of macrophages in suppressing the allergic effects has been described in a murine model of ovalbumin-induced allergic airway hyper-reactivity and DSS-induced colitis applying treatment with filarial cystatin[13]. Moreover, filarial cystatin reduces Respiratory Syncytial Virus (RSV)-inducing immunopathology by inducing high levels of IL-10 production by CD4<sup>+</sup> T cells in the airways and lungs of mice[41]. Cystatin from the nematode parasite *Ascaris lumbricoides* reduces inflammation in a mouse model of DSS-induced colitis by increasing the expression of IL-10, TGF $\beta$  with simultaneous reduction of IL-6 and TNF $\alpha$  [42]. Collectively, the studies described here indicate that cystatins from various nematodes not only function as immunomodulators in the infection setting but also have potential as therapeutic agents for inflammatory diseases. Our results highlight the unique role of cystatin to induce IL-10 in human monocytes and macrophages stimulated with microfilarial products. Also cystatin-dependent expression of PD-L1 and PD-L2 in human monocytes suggests that cystatin from microfilariae is partially involved in the immune regulation process and thus could contribute to asymptomatic infection through PD-L1:PD-1 dependent suppression of CD4<sup>+</sup> T cell function[26,43–45]. Since circulating microfilariae in asymptotically infected individuals will readily come into contact with blood monocytes, the modulation of this particular immune cell is vital for the parasite's immune evasion strategy.

Similar to our previously published results on Mf [26], stimulation of monocytes and macrophages with whole Mf resulted in the production of significantly higher levels of IL-8 compared to unstimulated controls. The significant elevation of IL-8 by Mf-stimulated human monocytes and macrophages, however, was not dependent on cystatin (Fig 3C and 3D). Thus, the effect by Mf could be attributed to the presence of other molecules in Mf that are also able to induce IL-8. So far there are only a few studies revealing the potential of filarial products to induce IL-8, such as lipoproteins from *Wolbachia*, an obligate intracellular symbiont of *Brugia malayi*, which was shown to induce IL-8 in a TLR2/6-dependent manner[46]. While infective L3 larvae are not able to induce IL-8 in dendritic cells (DC) [47] or epithelial cells[48], it is expressed in a co-culture of DC and keratinocytes[49]. In contrast to infective larvae, live Mf and excretory/secretory products of female worms can induce the expression of IL-8 in human DC[50] and monocytes[51]. Previously, and also in the present study, we could show significant elevation of IL-6 by Mf in monocytes only, but not in macrophages. Although we see an increased production of IL-6 in Mf-treated monocytes (Fig 3C), IL-6 was not significantly reduced in monocytes that were treated with  $\Delta$ Mf. This result suggests that other molecules from whole Mf lysate were responsible for the elevated levels of IL-6 by monocytes.

Furthermore, we studied the immune signalling events exploited by rBmCPI-2 leading to the production of IL-10 and IL-8 in human monocytes and macrophages. It has previously

been demonstrated that ES-62, an excretory-secretory product of the rodent filarial nematode *Acanthocheilonema viteae* induces tyrosine phosphorylation of glycoproteins in murine macrophages[52] and modulates the activation of MAP kinases (ERK, p38 and JNK), thereby regulating cytokine production[53,54]. A study from our own group revealed that IL-10 production in murine macrophages triggered by cystatin from *A. viteae* was tyrosine kinase sensitive, and depended on activation of both MAPK[35]. In addition, LPS-stimulated microglia of mice *in vitro* showed that increased IL-10 mRNA expressions by filarial cystatin was dependent on ERK signalling[55]. In agreement with this data, our results indicate a significant reduction of rBmCPI-2- or Mf-induced IL-10 in macrophages, suggesting roles for ERK and p38 in these cells. On the other hand, IL-10 induction by rBmCPI-2 or Mf in human monocytes was independent of ERK signalling, indicating a differential induction of IL-10 in this myeloid cell population. This suggests that filarial products are able to differentially exploit and/or trigger the signalling events in human monocytes and macrophages, leading to induction of IL-10. Moreover, while IL-10 was independent of the ERK pathway in human monocytes only, IL-8 production was dependent on both MAPK kinases (ERK and p38) in monocytes (Figs 5A and 6A) and macrophages (Figs 5B and 6B).

This is the first study addressing the contribution of filarial cystatin within microfilariae, which circulate in the blood, to induce IL-10 in monocytes. Overall, our results show that IL-10 and IL-8 induction by human filarial cystatin in human monocytes and macrophages is dependent on p38, while only IL-10 induction in macrophages is additionally dependent of ERK signalling. Moreover, Mf-induced IL-10 production, but not IL-8 or IL-6 induction in human monocytes and macrophages is dependent on cystatin, suggesting that while cystatin is sufficient and necessary for IL-10 induction, factors other than cystatin are responsible for IL-8 and IL-6.

## Supporting information

**S1 Fig. Induction of IL-10, but not IL-8 by microfilarial lysate (Mf) in human monocytes is independent of ERK.** Human monocytes (A) and macrophages (B) were left unstimulated or stimulated with 20 µg/ml of Mf for 24 h with various concentrations of MEK1/2 inhibitor. Cytokines in the supernatants were detected by ELISA. Data are from 7–9 donors and shown as mean ±SEM. Statistical analysis was done using the Wilcoxon matched-pairs signed rank test. Values statistically different from the unstimulated control are depicted as <sup>§</sup> p<0.05 and <sup>§§</sup> p<0.01. Values statistically different from Mf-stimulated monocytes are depicted as \* p<0.05 and \*\*p<0.01. (TIF)

**S2 Fig. Induction of IL-10 and IL-8 by Mf in human monocytes and macrophages is dependent on p38.** Human monocytes (A) and macrophages (B) were left unstimulated or stimulated with 20 µg/ml of Mf for 24 h with various concentrations of p38 inhibitor. Cytokines in the supernatants were detected by ELISA. Data are from 7–9 donors and shown as mean ±SEM. Statistical analysis was done using the Wilcoxon matched-pairs signed rank test. Values statistically different from the unstimulated control are depicted as <sup>§</sup> p<0.05 and <sup>§§</sup> p<0.01. Values statistically different from Mf-stimulated monocytes are depicted as \* p<0.05 and \*\*p<0.01. (TIF)

## Acknowledgments

We thank the NIAID/NIH Filariasis Research Reagent Resource Center (FR3) in Athens, Georgia for the kind donation of *Brugia malayi* microfilariae. We thank Bettina Sonnenburg

for her excellent technical assistance and Ankur Midha and Ivet Yordanova for critically reviewing the manuscript.

## Author Contributions

**Conceptualization:** Susanne Hartmann, Svenja Steinfelder.

**Data curation:** Gopinath Venugopal, Susanne Hartmann.

**Formal analysis:** Gopinath Venugopal, Svenja Steinfelder.

**Funding acquisition:** Susanne Hartmann.

**Investigation:** Gopinath Venugopal, Marion Mueller, Svenja Steinfelder.

**Methodology:** Gopinath Venugopal, Marion Mueller, Svenja Steinfelder.

**Project administration:** Susanne Hartmann, Svenja Steinfelder.

**Resources:** Susanne Hartmann.

**Software:** Susanne Hartmann.

**Supervision:** Susanne Hartmann, Svenja Steinfelder.

**Validation:** Susanne Hartmann, Svenja Steinfelder.

**Visualization:** Gopinath Venugopal, Svenja Steinfelder.

**Writing ± original draft:** Gopinath Venugopal.

**Writing ± review & editing:** Susanne Hartmann, Svenja Steinfelder.

## References

1. McSorley HJ, Maizels RM. Helminth infections and host immune regulation. *Clin Microbiol Rev.* 2012; 25: 585±608. <https://doi.org/10.1128/CMR.05040-11> PMID: 23034321
2. Hotez PJ, Brindley PJ, Bethony JM, King CH, Pearce EJ, Jacobson J. Helminth infections: the great neglected tropical diseases. *J Clin Invest.* 2008; 118: 1311±1321. <https://doi.org/10.1172/JCI34261> PMID: 18382743
3. Anthony RM, Rutitzky LI, Urban JF Jr., Stadecker MJ, Gause WC. Protective immune mechanisms in helminth infection. *Nat Rev Immunol.* 2007; 7: 975±987. <https://doi.org/10.1038/nri2199> PMID: 18007680
4. Allen JE, Maizels RM. Diversity and dialogue in immunity to helminths. *Nat Rev Immunol.* Nature Publishing Group; 2011; 11: 375±88. <https://doi.org/10.1038/nri2992> PMID: 21610741
5. Johnston KL, Ford L, Taylor MJ. Overcoming the challenges of drug discovery for neglected tropical diseases: the A.WOL experience. *J Biomol Screen.* 2014; 19: 335±343. <https://doi.org/10.1177/1087057113511270> PMID: 24241712
6. World Health Organization. Global Health Estimates 2015: Deaths by Cause, Age, Sex, by Country and by Region, 2000±2015 [Internet]. Global Health Estimates: Key figures and tables. 2016. Available: [http://www.who.int/healthinfo/global\\_burden\\_disease/estimates/en/index1.html](http://www.who.int/healthinfo/global_burden_disease/estimates/en/index1.html)
7. Fenwick A. The global burden of neglected tropical diseases. *Public Health.* Elsevier Ltd; 2012; 126: 233±236. <https://doi.org/10.1016/j.puhe.2011.11.015> PMID: 22325616
8. Babu S, Nutman TB. Immunology of lymphatic filariasis. *Parasite Immunol.* 2014; 36: 338±346. <https://doi.org/10.1111/pim.12081> PMID: 24134686
9. Hartmann S, Kyewski B, Sonnenburg B, Lucius R. A filarial cysteine protease inhibitor down-regulates T cell proliferation and enhances interleukin-10 production. *Eur J Immunol.* 1997; 27: 2253±2260. <https://doi.org/10.1002/eji.1830270920> PMID: 9341767
10. Manoury B, Gregory WF, Maizels RM, Watts C. Bm-CPI-2, a cystatin homolog secreted by the filarial parasite *Brugia malayi*, inhibits class II MHC-restricted antigen processing. *Curr Biol.* 2001; 11: 447±451. [https://doi.org/10.1016/S0960-9822\(01\)00118-X](https://doi.org/10.1016/S0960-9822(01)00118-X) PMID: 11301256

11. Dainichi T, Maekawa Y, Ishii K, Zhang T, Fawzy Nashed B, Sakai T, et al. Nippocystatin, a cysteine protease inhibitor from *Nippostrongylus brasiliensis*, inhibits antigen processing and modulates antigen-specific immune response. *Infect Immun*. 2001; 69: 7380±7386. <https://doi.org/10.1128/IAI.69.12.7380-7386.2001> PMID: 11705911
12. Schenemeyer A, Lucius R, Sonnenburg B, Brattig N, Sabat R, Schilling K, et al. Modulation of human T cell responses and macrophage functions by onchocystatin, a secreted protein of the filarial nematode *Onchocerca volvulus*. *J Immunol*. 2001; 167: 3207±15. doi:11544307 PMID: 11544307
13. Schnoeller C, Rausch S, Pillai S, Avagyan A, Wittig BM, Loddenkemper C, et al. A helminth immunomodulator reduces allergic and inflammatory responses by induction of IL-10-producing macrophages. *J Immunol*. 2008; 180: 4265±72. doi:180/6/4265 [pii] PMID: 18322239
14. Pfaff AW, Schulz-Key H, Soboslay PT, Taylor DW, MacLennan K, Hoffmann WH. Litomosoides sigmondontis cystatin acts as an immunomodulator during experimental filariasis. *Int J Parasitol*. 2002; 32: 171±178. [https://doi.org/10.1016/S0020-7519\(01\)00350-2](https://doi.org/10.1016/S0020-7519(01)00350-2) PMID: 11812494
15. Gomez-Escobar N, Gregory WF, Maizels RM. Identification of tgh-2, a filarial nematode homolog of *Caenorhabditis elegans* daf-7 and human transforming growth factor??, expressed in microfilarial and adult stages of *Brugia malayi*. *Infect Immun*. 2000; 68: 6402±6410. <https://doi.org/10.1128/IAI.68.11.6402±6410.2000> PMID: 11035752
16. Gomez-Escobar N, Lewis E, Maizels RM. A Novel Member of the Transforming Growth Factor-β (TGF-β) Superfamily from the Filarial Nematode *Brugia malayi* and *B. pahangi*. *Exp Parasitol*. 1998; 88: 200±209. <https://doi.org/10.1006/expr.1998.4248> PMID: 9562423
17. Grainger JR, Smith KA, Hewitson JP, McSorley HJ, Harcus Y, Filbey KJ, et al. Helminth secretions induce de novo T cell Foxp3 expression and regulatory function through the TGF-β pathway. *J Exp Med*. 2010; 207: 2331±41. <https://doi.org/10.1084/jem.20101074> PMID: 20876311
18. Zang X. Homologues of Human Macrophage Migration Inhibitory Factor from a Parasitic Nematode. GENE CLONING, PROTEIN ACTIVITY, AND CRYSTAL STRUCTURE. *J Biol Chem*. 2002; 277: 44261±44267. <https://doi.org/10.1074/jbc.M204655200> PMID: 12221083
19. Hewitson JP, Grainger JR, Maizels RM. Helminth immunoregulation: The role of parasite secreted proteins in modulating host immunity. *Mol Biochem Parasitol*. 2009; 167: 1±11. <https://doi.org/10.1016/j.molbiopara.2009.04.008> PMID: 19406170
20. Hartmann S, Lucius R. Modulation of host immune responses by nematode cystatins. *Int J Parasitol*. 2003; 33: 1291±1302. [https://doi.org/10.1016/S0020-7519\(03\)00163-2](https://doi.org/10.1016/S0020-7519(03)00163-2) PMID: 13678644
21. Grzonka Z, Jankowska E, Kasprzykowski F, Kasprzykowska R, Rankiewicz L, Wiczek W, et al. Structural studies of cysteine proteases and their inhibitors. *Acta Biochim Pol*. 2001; 48: 1±20. [https://doi.org/10.1016/0300-9629\(90\)90008-G](https://doi.org/10.1016/0300-9629(90)90008-G) PMID: 11440158
22. Rzychon M, Chmiel D, Stec-Niemczyk J. Modes of inhibition of cysteine proteases. *Acta Biochim Pol*. 2004; 51: 861±873. doi:045104861 <https://doi.org/045104861> PMID: 15625558
23. Sabotič J, Kos J. Microbial and fungal protease inhibitors: Current and potential applications. *Appl Microbiol Biotechnol*. 2012; 93: 1351±1375. <https://doi.org/10.1007/s00253-011-3834-x> PMID: 22218770
24. Kopitar-Jerala N. The role of cysteine proteinases and their inhibitors in the host-pathogen cross talk. *Curr Protein Pept Sci*. 2012; 13: 767±75. <https://doi.org/10.2174/138920312804871102> PMID: 23305363
25. Hartmann S, Schenemeyer A, Sonnenburg B, Vray B, Lucius R. Cystatins of filarial nematodes up-regulate the nitric oxide production of interferon-γ-activated murine macrophages. 2002; 253±262.
26. O'Regan NL, Steinfeldt S, Venugopal G, Rao GB, Lucius R, Srikantam A, et al. *Brugia malayi* Microfilariae Induce a Regulatory Monocyte/Macrophage Phenotype That Suppresses Innate and Adaptive Immune Responses. *PLoS Negl Trop Dis*. 2014; 8. <https://doi.org/10.1371/journal.pntd.0003206> PMID: 25275395
27. Babu S, Blauvelt CP, Kumaraswami V, Nutman TB. Diminished expression and function of TLR in lymphatic filariasis: a novel mechanism of immune dysregulation. *J Immunol*. 2005; 175: 1170±1176. doi:175/2/1170 [pii] PMID: 16002719
28. Semnani RT, Keiser PB, Coulibaly YI, Keita F, Diallo AA, Traore D, et al. Filaria-induced monocyte dysfunction and its reversal following treatment. *Infect Immun*. 2006; 74: 4409±4417. <https://doi.org/10.1128/IAI.01106-05> PMID: 16861626
29. Babu S, Kumaraswami V, Nutman TB. Alternatively activated and immunoregulatory monocytes in human filarial infections. *J Infect Dis*. 2009; 199: 1827±37. <https://doi.org/10.1086/599090> PMID: 19456233
30. Asadullah K, Sterry W, Volk HD. Interleukin-10 Therapy: Review of a New Approach. *Med Immunol*. 2003; 55: 241±269. <https://doi.org/10.1124/pr.55.2.4.241>

31. Satapathy AK, Sartono E, Sahoo PK, Dentener MA, Michael E, Yazdanbakhsh M, et al. Human bancroftian filariasis: immunological markers of morbidity and infection. *Microbes Infect.* 2006; 8: 2414±2423. <https://doi.org/10.1016/j.micinf.2006.05.003> PMID: 16839794
32. Babu S, Blauvelt CP, Kumaraswami V, Nutman TB. Chemokine receptors of T cells and of B cells in lymphatic filarial infection: a role for CCR9 in pathogenesis. *J Infect Dis.* 2005; 191: 1018±1026. <https://doi.org/10.1086/427658> PMID: 15717282
33. Gregory WF, Maizels RM. Cystatins from filarial parasites: Evolution, adaptation and function in the host-parasite relationship. *Int J Biochem Cell Biol.* 2008; 40: 1389±1398. <https://doi.org/10.1016/j.biocel.2007.11.012> PMID: 18249028
34. Bennuru S, Semnani R, Meng Z, Ribeiro JMC, Veenstra TD, Nutman TB. Brugia malayi excreted/secreted proteins at the host/parasite interface: Stage- and gender-specific proteomic profiling. *PLoS Negl Trop Dis.* 2009; 3. <https://doi.org/10.1371/journal.pntd.0000410> PMID: 19352421
35. Klotz C, Ziegler T, Figueiredo AS, Rausch S, Hepworth MR, Obsivac N, et al. A helminth immunomodulator exploits host signaling events to regulate cytokine production in macrophages. *PLoS Pathog.* 2011; 7. <https://doi.org/10.1371/journal.ppat.1001248> PMID: 21253577
36. Maizels RM, Yazdanbakhsh M. Immune regulation by helminth parasites: cellular and molecular mechanisms. *Nat Rev Immunol.* 2003; 3: 733±744. <https://doi.org/10.1038/nri1183> PMID: 12949497
37. Specht S, Taylor MD, Hoeve MA, Allen JE, Lang R, Hoerauf A. Over expression of IL-10 by macrophages overcomes resistance to murine filariasis. *Exp Parasitol.* Elsevier Inc.; 2012; 132: 90±96. <https://doi.org/10.1016/j.exppara.2011.09.003> PMID: 21959021
38. Figueiredo CA, Barreto ML, Rodrigues LC, Cooper PJ, Silva NB, Amorim LD, et al. Chronic intestinal helminth infections are associated with immune hyporesponsiveness and induction of a regulatory network. *Infect Immun.* 2010; 78: 3160±3167. <https://doi.org/10.1128/IAI.01228-09> PMID: 20404082
39. Araujo MIAS, Hoppe B, Medeiros M, Alca L, Cec m, Schriefer A, et al. Impaired T Helper 2 Response to Aeroallergen in Helminth-Infected Patients with Asthma. *J Infect Dis.* 2004; 190: 1797±1803. <https://doi.org/10.1086/425017> PMID: 15499536
40. Turner JD, Jackson JA, Faulkner H, Behnke J, Else KJ, Kamgno J, et al. Intensity of intestinal infection with multiple worm species is related to regulatory cytokine output and immune hyporesponsiveness. *J Infect Dis.* 2008; 197: 1204±1212. <https://doi.org/10.1086/586717> PMID: 18462167
41. Schuijs MJ, Hartmann S, Selkirk ME, Roberts LB, Openshaw PJM, Schnoeller C. The Helminth-derived immunomodulator AvCystatin reduces virus enhanced inflammation by induction of regulatory IL-10+ T cells. *PLoS One.* 2016; 11: 1±14. <https://doi.org/10.1371/journal.pone.0161885> PMID: 27560829
42. Coronado S, Barrios L, Zakzuk J, Regino R, Ahumada V, Franco L, et al. A recombinant cystatin from *Ascaris lumbricoides* attenuates inflammation of DSS-induced colitis. *Parasite Immunol.* 2017; e12425. <https://doi.org/10.1111/pim.12425> PMID: 28295446
43. Freeman GJ, Long AJ, Iwai Y, Bourque K, Chernova T, Nishimura H, et al. Engagement of the Pd-1 Immunoinhibitory Receptor by a Novel B7 Family Member Leads to Negative Regulation of Lymphocyte Activation. *J Exp Med.* 2000; 192: 1027±1034. <https://doi.org/10.1084/jem.192.7.1027> PMID: 11015443
44. Butte MJ, Keir ME, Phamduy TB, Sharpe AH, Freeman GJ. Programmed Death-1 Ligand 1 Interacts Specifically with the B7-1 Costimulatory Molecule to Inhibit T Cell Responses. *Immunity.* 2007; 27: 111±122. <https://doi.org/10.1016/j.immuni.2007.05.016> PMID: 17629517
45. Wilcox RA, Feldman AL, Wada DA, Yang ZZ, Comfere NI, Dong H, et al. B7-H1 (PD-L1, CD274) suppresses host immunity in T-cell lymphoproliferative disorders. *Blood.* 2009; 114: 2149±2158. <https://doi.org/10.1182/blood-2009-04-216671> PMID: 19597183
46. Turner JD, Langley RS, Johnston KL, Gentil K, Ford L, Wu B, et al. Wolbachia lipoprotein stimulates innate and adaptive immunity through toll-like receptors 2 and 6 to induce disease manifestations of filariasis. *J Biol Chem.* 2009; 284: 22364±22378. <https://doi.org/10.1074/jbc.M901528200> PMID: 19458089
47. Semnani RT, Law M, Kubofcik J, Nutman TB. Filaria-Induced Immune Evasion: Suppression by the Infective Stage of Brugia malayi at the Earliest Host-Parasite Interface. *J Immunol.* 2004; 172: 6229±6238. <https://doi.org/10.4049/jimmunol.172.10.6229> PMID: 15128811
48. Schroeder JH, Simbi BH, Ford L, Cole SR, Taylor MJ, Lawson C, et al. Live Brugia malayi Microfilariae Inhibit Transendothelial Migration of Neutrophils and Monocytes. *PLoS Negl Trop Dis.* 2012; 6. <https://doi.org/10.1371/journal.pntd.0001914> PMID: 23209856
49. Boyd A, Bennuru S, Wang Y, Sanprasert V, Law M, Chaussabel D, et al. Quiescent Innate Response to Infective Filariae by Human Langerhans Cells Suggests a Strategy of Immune Evasion. *Infect Immun.* 2013; 81: 1420±1429. <https://doi.org/10.1128/IAI.01301-12> PMID: 23429540

50. Semnani RT, Liu AY, Sabzevari H, Kubofcik J, Zhou J, Gilden JK, et al. *Brugia malayi* microfilariae induce cell death in human dendritic cells, inhibit their ability to make IL-12 and IL-10, and reduce their capacity to activate CD4+ T cells. *J Immunol*. 2003; 171: 1950±60. <https://doi.org/10.4049/jimmunol.171.4.1950> PMID: 12902498
51. Weinkopff T, Mackenzie C, Eversole R, Lammie PJ. Filarial Excretory-Secretory Products Induce Human Monocytes to Produce Lymphangiogenic Mediators. *PLoS Negl Trop Dis*. 2014; 8. <https://doi.org/10.1371/journal.pntd.0002893> PMID: 25010672
52. Goodridge HS, Deehan MR, Harnett W, Harnett MM. Subversion of immunological signalling by a filarial nematode phosphorylcholine-containing secreted product. *Cell Signal*. 2005; 17: 11±16. <https://doi.org/10.1016/j.cellsig.2004.05.014> PMID: 15451020
53. Goodridge HS, Harnett W, Liew FY, Harnett MM. Differential regulation of interleukin-12 p40 and p35 induction via Erk mitogen-activated protein kinase-dependent and -independent mechanisms and the implications for bioactive IL-12 and IL-23 responses. *Immunology*. 2003; 109: 415±425. <https://doi.org/10.1046/j.1365-2567.2003.01689.x> PMID: 12807488
54. Harnett W, Goodridge HS, Harnett MM. Subversion of immune cell signal transduction pathways by the secreted filarial nematode product, ES-62. *Parasitology*. 2005; 130 Suppl: S63±8. <https://doi.org/10.1017/S0031182005008164> PMID: 16281993
55. Behrendt P, Arnold P, Brueck M, Rickert U, Lucius R, Hartmann S, et al. A Helminth Protease Inhibitor Modulates the Lipopolysaccharide-Induced Proinflammatory Phenotype of Microglia in vitro. *Neuroimmunomodulation*. 2016; 23: 109±121. <https://doi.org/10.1159/000444756> PMID: 27088850