

IL-4

Interleukin 4, B-cell stimulatory factor murine, recombinant, *E. coli*

Cat. No.	Amount
PR-464	10 µg

For *in vitro* use only
Quality guaranteed for 12 months
Store at -20°C

Avoid freeze / thaw cycles

Form

Lyophilized.

Solubility

It is recommended to reconstitute the lyophilized IL-4 in bidest H₂O not less than 100 µg/ml, which can then be further diluted to other aqueous solutions. For long term storage it is recommended to add a carrier protein (0.1% HSA or BSA).

Activity

ED₅₀: < 2 ng/ml, determined by the dose-dependent proliferation of murine HT-2 cells.

Endotoxin

Less than 0.1 ng/µg (IEU/µg) of IL-4.

Molecular Weight

13.5 kDa

Purity

≥ 95% by SDS-PAGE and RP-HPLC

Description

The anti-inflammatory cytokine Interleukin-4 is involved in the regulation of inflammatory responses by promoting the differentiation of naïve T-cells to T helper type 2 (Th2) cells. A key anti-inflammatory action of IL-4 results from its ability to inhibit the release of pro-inflammatory cytokines by innate immune cells and to upregulate the synthesis of IL-1 receptor antagonist.

IL-4 has been demonstrated to be a potent cofactor for B and T lymphocyte proliferation and differentiation. Recombinant murine IL-4 produced in *E. coli* is a single, non-glycosylated form of murine IL-4 polypeptide chain containing 120 amino acids and having a molecular mass of 13.5 kDa. IL-4 is purified by proprietary chromatographic techniques.

Selected References:

- Wurster *et al.* (2002) Interleukin-4-mediated protection of primary B cells from apoptosis through Stat6-dependent up-regulation of BclxL. *J. Biol. Chem.* **277**:27169.
- Nolan *et al.* (2004) Role of interleukin-4 in regulation of age-related inflammatory changes in the hippocampus. *J. Biol. Chem.* **280**:9354.
- McCurdy *et al.* (2004) Modified vaccinia virus Ankara immunization protects against lethal challenge with recombinant vaccinia virus expressing murine interleukin-4. *J. Virol.* **78**:12471.
- Yatsenko *et al.* (2004) Alternative splicing of murine interleukin-4 mRNA. *Bull. Exp. Biol. Med.* **137**:179.
- Walz *et al.* (2002) A murine interleukin-4-Ig fusion protein regulates the expression of Th1- and Th2-specific cytokines in the pancreas of NOD mice. *Horm. Metab. Res.* **34**:561.
- Ghiasi *et al.* (2001) Recombinant herpes simplex virus type 1 expressing murine interleukin-4 is less virulent than wild-type virus in mice. *J. Virol.* **75**:9029.
- Yoshikawa *et al.* (2000) Immune gene therapy of experimental mouse brain tumor with adenovirus-mediated gene transfer of murine interleukin-4. *Cancer Immunol. Immunother.* **49**:23.
- Missol-Kolka *et al.* (1998) Combined therapy of B16(F10) murine melanoma using *E. coli* cytosine deaminase gene and murine interleukin-4 gene. *Neoplasma.* **45**:305.