# Bioluminescent *Leishmania* expressing luciferase for rapid and high throughput screening of drugs acting on amastigote-harbouring macrophages and for quantitative real-time monitoring of parasitism features in living mice

# Thierry Lang,\* Sophie Goyard, Mai Lebastard and Geneviève Milon

Unité d'Immunophysiologie et Parasitisme Intracellulaire, Département de Parasitologie, Institut Pasteur, rue 25 du Dr Roux, 75724 Paris cedex 15, France.

## **Summary**

In this study, we have established conditions for generating Leishmania amazonensis recombinants stably expressing the firefly luciferase gene. These parasites produced significant bioluminescent signals for both in vitro studies and the development of an in vivo model, allowing the course of the parasitism to be readily monitored in real time in the living animals such as laboratory mice. First, a model was established, using parasite-infected mouse macrophages for rapidly determining the activity of drugs against intracellular amastigotes. Results indicated that recombinant Leishmania can be reliably and confidently used to monitor compounds acting on intracellular amastigote-harbouring macrophages. Secondly, temporal analyses were performed following inoculation of metacyclic promastigotes into the ear dermis of BALB/c mice and the bioluminescent light transmitted through the tissue was imaged externally using a charge coupled device (CCD) camera. Bioluminescent signals, measured at the inoculation site and in the draining lymph node of mice containing these parasites correlated well with the more classical quantification of parasites. These assays prove that the real-time bioluminescent assay is not only sensitive but also more rapid than culture-base techniques allowing to monitor parasite-load before any clinical signs of leishmaniasis are detectable. In short, this luciferase imaging study is useful to monitor the efficacy of anti-leishmanial drugs on live cell culture and to trace leishmanial infection in animal models.

Received 9 July, 2004; revised 9 September, 2004; accepted 10 September, 2004. \*For correspondence. E-mail tlang@pasteur.fr; Tel. (+33) 1 45 68 86 73; Fax (+33) 1 40 61 33 32.

#### Introduction

Leishmania spp. are intracellular protozoan parasites transmitted from the sandfly during a blood meal to a variety of mammalian hosts. They cause a large spectrum of diseases affecting the skin mucous membranes and viscera. These dimorphic parasites exist as extracellular flagellated promastigotes in the insect vector and as obligatory intracellular amastigotes in the mammalian host. Experimental models, such as in vitro culture of mouse- or human-derived macrophages or laboratory mice, are largely used to study immuno-biology of these parasites and to screen the efficacy of newly developed drugs as well as vaccines (Blackwell, 1996; Launois et al., 1997; Handman, 2001). Simple, reliable and rapid test allowing the determination of drug activity at the mammalian stage are still needed for the development of new drugs and require detection and quantification of the Leishmania amastigote burdens in different host cells. This is routinely performed by Giemsa staining (Berman, 1984), culture-based techniques (Titus et al., 1985; Buffet et al., 1995), classic polymerase chain reaction (PCR)based methods (Katakura et al., 1998; Noyes et al., 1998) or real-time quantitative PCR (Nicolas et al., 2002). However these techniques are cumbersome, laborious and rendered difficult the screening of several drugs at a time. Furthermore, while some of these methods are useful for in vitro studies, they have not proved ideal for the understanding of biological processes as they occur in living animals.

To better investigate parasite expansion, both *ex vivo* and *in vivo* in living cells, rapid, non-destructive, real-time monitoring methods that are adaptable to the experimental animal model are needed. To this end, reporter genes offer a method of labelling invasive microorganisms that is innocuous and allows the sensitive detection of only live, metabolically active cells by biophotonic imaging. The firefly luciferase represents one of the most efficient biological reporters. It catalyses the reaction of the substrate luciferin with adenosine triphosphate (ATP) to generate photons. Monitoring pathogenic processes in a living animal was first demonstrated with strains of Sal-

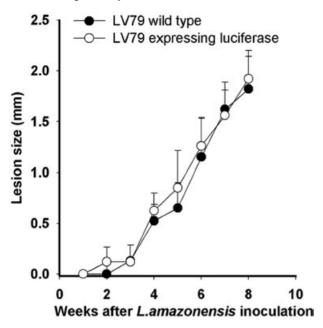


Fig. 1. Course of lesion development in BALB/c mice intradermally inoculated with 104 metacyclic promastigotes of L. amazonensis-wildtype strain or -expressing luciferase.

monella (Contag et al., 1995). This non-invasive, rapid, real-time monitoring has been applied to various bacterial infection models and has been shown to have significant advantages other conventional methods for studying microorganism dissemination and treatment in animals (Loimaranta et al., 1998; Francis et al., 2000; 2001; Rocchetta et al., 2001; Edinger et al., 2003) or monitoring activities of bacteria within biofilm in a mouse model (Kadurugamuwa et al., 2003). The use of luciferase gene has also improved the in vitro screening of antimicrobial agents acting on microorganisms developing within host cells (Deb et al., 2000). The same strategy has been of great interest while testing new anti-leishmanial agents under both axenic and promastigote conditions (Luque-Ortega et al., 2001; Sereno et al., 2001).

In this study, we have established conditions for transfection of Leishmania amazonensis recombinants stably expressing the firefly luciferase gene. This luciferase expressing parasite was further analysed in two different systems: (i) in vitro drug screening on the clinically relevant stage of the parasites, i.e. amastigote-loaded mouse bone marrow derived macrophages, and (ii) the monitoring of the parasitic process in living animals. Thus, the light-emitting metacyclic promastigotes were injected into the ear dermis of BALB/c mice and the bioluminescent light transmitted through the tissue was imaged externally using a charge coupled device (CCD) camera at different time points after parasite delivery.

#### Results

Onset and development of lesions in BALB/c mice infected with bioluminescent L. amazonensis

Luciferase expressing-metacyclic promastigotes were injected into the ear dermis of BALB/c mice. The development of lesions was assessed by measuring the thickness of the ear over a period of 8 weeks (Fig. 1) and compared with the lesions induced by the inoculation of the same dose of wild-type promastigotes. In both cases, the onset of cutaneous clinical signs was detectable around week 3 and exhibiting ear lesions from week 7, which then steadily increased in size during the period examined (Fig. 1). This experiment clearly indicated that both control and luciferase-expressing parasites are driving the same clinical processes in BALB/c mice.

Bioluminescence assay on parasite populations at various developmental stages

To assess the feasibility of using bioluminescence as a quantitative indicator of the viability and the multiplication of parasites, studies were performed to compare bioluminescence to the number of viable parasites. Wild-type strain of Leishmania was used as control. After washing the cells, serial dilutions of parasites  $(2.5 \times 10^6 - 1.2 \times 10^3)$ were made to measure luciferase activity (Fig. 2).

The maximum level of bioluminescence was reached 5 min after substrate addition, followed by slow decay

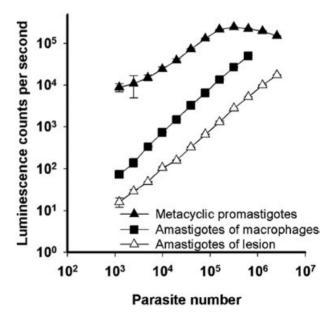


Fig. 2. Luminescence assay on luciferase-transfected L. amazonensis. Results for metacyclic promastigotes, amastigotes from BALB/c lesion or from infected macrophages are shown. The bioluminescent reaction was carried out as described in Experimental procedures with the Luciferase Reporter Gene Assay and the light emission was measured with the luminometer.

during 1 h. No bioluminescence (<10 cps) was observed in the wild-type strain at any the parasite stage (not shown), indicating that bioluminescence is specific to the luciferase-expressing parasites.

During log phase growth the luminescence signal was proportional to the number of bioluminescent parasites (data not shown). Once the cultures reached the stationary phase the bioluminescence decreased. This data suggests that a decrease in growth correlated with a decrease in metabolic activity of the population. When assayed with metacyclic promastigotes, luminescence  $(8.8 \times 10^3)$ counts per second) was detected for  $1.2 \times 10^3$  parasites. proportional to the cell number and reached  $1.3 \times 10^5$  cps for  $7.8 \times 10^4$  parasites (Fig. 2). We next investigated the bioluminescence of lesion-derived parasites and compared it with amastigotes purified from infected macrophages. Amastigotes were purified either from the L. amazonensis-infected footpads or from the in vitro Leishmania-harbouring macrophages. The luminescence obtained with lesion-derived amastigotes was 3-6 times lower than with amastigotes of macrophages and 10-50 times lower than with metacyclic promastigotes. The minimum level (47 cps for 1.2 × 10<sup>3</sup> amastigotes) was 10 times higher than the level measured with the wild-type strain. These results indicate that all the parasites recovered from lesions were luminescent confirming the stability of luciferase-transfected parasites in this type of lesion.

Validation of the luminescence assay using Leishmaniainfected macrophages: monitoring with drugs known to be active on Leishmania spp.

Bioluminescence of Leishmania was measured in infected macrophages and inhibition was monitored in presence of anti-leishmanial drugs. Macrophages were infected with stationary-phase culture of transfected-Leishmania at an 1:8 multiplicity of infection in 24-well plate. The fate of Leishmania was then followed over a period of 48 h by measuring luminescence as well as the percentage of infected macrophages. The transfectants showed characteristics of the parental cells. They were infective for the macrophages and the differentiation into amastigotes with typical parasitophorous vacuole seen for *L. amazonensis* (Antoine et al., 1991) was observed. The inhibition of growth by the anti-leishmanial compound amphotericin B was demonstrated. Over a period of 48 h, the percentage of infected macrophages at the time that the drug was added was high (>85%) and remained constant when no drug was added to the medium (Fig. 3A). However, after the treatment of macrophages with amphotericin, the percentage of infected cells and intensity of bioluminescence decreased proportionally to the drug concentration (Fig. 3A and B).

The high intensity and stability of bioluminescence of the luciferase-expressing L. amazonensis amastigotes

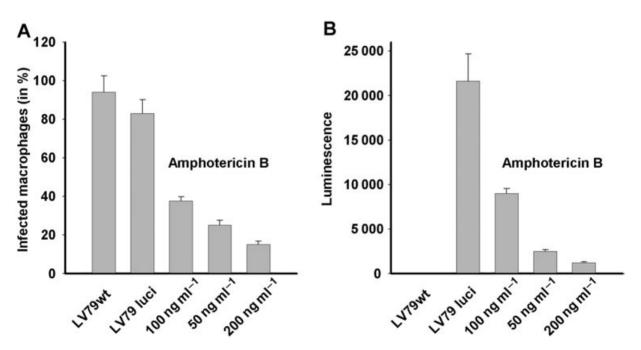


Fig. 3. Percentage (A) and bioluminescence intensity (B) of Leishmania-infected macrophages treated or not with amphotericin. Bone marrowderived macrophages were prepared from BALB/c mice and infected with luciferase-transfected L. amazonensis promastigotes (eight parasites per host cell). Wild-type strain was used as negative control. Two hours later extracellular parasites were removed and after 24 h amphotericin B was added (200, 100, 50 ng ml<sup>-1</sup>). The luciferase activity and the percentage of infected was measured 48 h after infection. The bioluminescent reaction in cells was triggered by the addition of the reagents for the determination of luciferase activity of the firefly, and the light emission was measured with a microtitration plate luminometer. The results are the mean of duplicates and representative of two experiments.

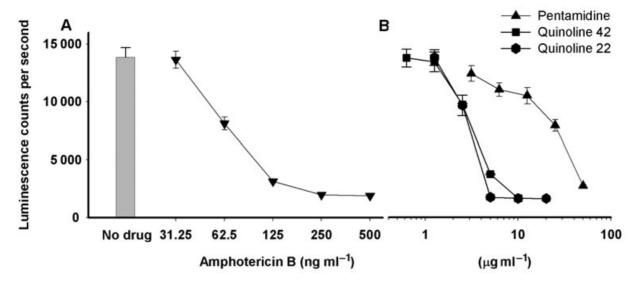


Fig. 4. Macrophages  $(4 \times 10^4)$  were infected in 96-well plate with recombinant *Leishmania* and treated with amphotericin (A) or pentamidine (B) or quinolines (B). The drugs were added as described in Fig. 3 and the luciferase activity of infected cells was measured at 48 h with the Luciferase Reporter Gene Assay and the light emission was measured with the luminometer. The results are the mean of duplicates and representative of three experiments.

allowed drug testing in microplates. The inhibition of growth by the anti-leishmanial compounds amphotericin B, pentamidine or quinolines was demonstrated (Fig. 4). In both cases the difference in the level of luminescence corresponded to a difference in the percentage of infected cells, observed microscopically (not shown). These findings show that luciferase-expressing *Leishmania* could be useful in a bioluminescence assay for drug screening in 96-well plates.

Ex vivo bioluminescence imaging of tissues from BALB/c mice inoculated with L. amazonensis in the footpad several months earlier

After inoculation of  $10^6$  parasites in the footpad, small nodular lesions were detected in the mice from week 3; they increased quickly in size and displayed ulceration and tissue necrosis after 12 weeks. The onset of the cutaneous lesions was associated with a large increase in the number of parasites in the infected footpad ( $>5 \times 10^8$  parasites). By contrast the number of parasites per draining lymph node remained low during the first weeks post-inoculation (about 100 parasites) and increased slowly in parallel with lesion progression until week 24 ( $10^4$  parasites). No viable parasites could be recovered from the spleens during the onset of lesions and less than 500 parasites per spleen were detected by week 24 post-inoculation.

Mice primarily infected 24 weeks before in the footpad with 10<sup>6</sup> parasites received luciferin and were sacrificed. The tissues were surgically removed and bioluminescence signal was analysed at the site of inoculation and its draining lymph node as well as distant tissues including

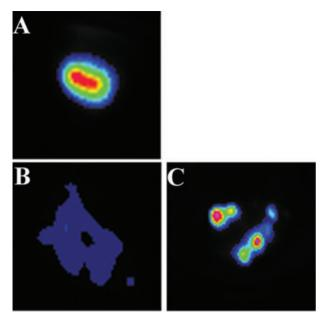
the spleen. Figure 5A illustrates the luminescent image of a representative lymph node; the bioluminescent signal was clearly detected and reached approximately  $3\times10^8$  photons per second (Ph/s). By contrast the intensity of the bioluminescent signal in the footpad appeared to be lower  $(4\times10^7$  Ph/s; Fig. 5B) than that in the lymph node which could be due to the lack of both oxygen and luciferin to the wounded tissue. This last point was confirmed by the fact that the footpad exhibited a higher luminescence activity  $(3\times10^8$  Ph/s) when luciferin was added a second time in liquid medium (Fig. 5C). In the spleen the luminescence was not detectable.

In vivo bioluminescence of L. amazonensis metacyclic promastigotes in living animals

We next investigated the parasite burden during the first 2 weeks after parasite inoculation by determining the number of parasites present in infected ears and expressing them as limiting dilution assay units (LDAU) (Fig. 6). One day after the injection of 10<sup>6</sup> parasites, viable parasites were detected in the ear (about 2% of the inoculum). The number of parasites then increased slowly during the first week after inoculation and then increased substantially reaching 10<sup>6</sup> parasites by day 14. Successful infection, defined as the development of lesion in ear dermis, was observed in 100% of the inoculated mice.

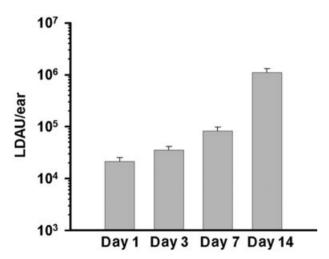
Figures 7 and 8 illustrate the real-time luminescent images from representative mice, infected with luciferase-transfected parasites over a 16 day time-course. Mice were infected with 10<sup>6</sup>, 10<sup>5</sup> or 10<sup>4</sup> parasites in the ear. Parasites produced a significant bioluminescent signal in

© 2004 Blackwell Publishing Ltd, Cellular Microbiology, 7, 383-392

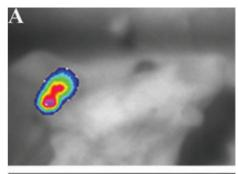


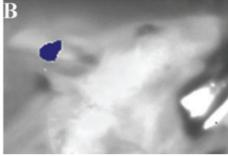
**Fig. 5.** Ex vivo imaging luminescence in the lymph node (A) and footpad (B, C) of BALB/c mice subcutaneously inoculated in the footpad with  $1 \times 10^6$  luciferase-transfected *Leishmania amazonensis*. Six months after infection, luciferin was intraperitoneally injected in mice. Mice were then sacrificed and the lymph node and the footpad were removed before analysing by luminescence. In C, luciferin was added a second time directly in the medium. The bioluminescence was monitored by detecting photon emission using a CCD camera. The bioluminescent signal is presented as a pseudo-colour image representing light intensities (red most and blue least intense).

mice allowing the progression of infection to be monitored non-invasively in the experimental model. The total photon emission from the infected site was quantified and shown in Fig. 8. No bioluminescence was observed in uninfected



**Fig. 6.** Monitoring of parasite load in BALB/c mice inoculated intradermally in ear with 10<sup>6</sup> metacyclic promastigotes of *Leishmania amazonensis*. At various time-points after inoculation, parasite loads were estimated in the ears by limiting dilution assays. Results are expressed in LDAU per ear. Histograms show the mean values + one standard deviation for three individual ears.



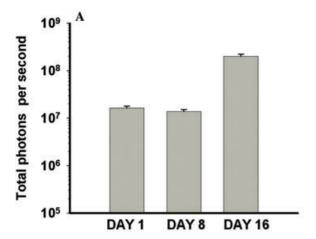


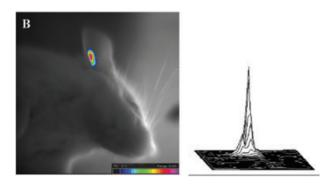
**Fig. 7.** *In vivo* luminescence imaging of *Leishmania* in a mouse model. Luciferase-transfected *Leishmania amazonensis* were inoculated at intradermal site with doses ranging from  $1\times10^6$  (A) to  $1\times10^5$  (B) and detection in the ear dermis was monitored by detecting photon emission at day 1 using a CCD camera. The bioluminescent signal is presented as a pseudo-colour image representing light intensities over the body surface area. Red represents the most intense light emission while blue correspond to the weakest signal. This pseudo-image is superimposed on the photography for orientation.

mice or mice infected with the wild-type strain. Following inoculation of transfected parasites, the bioluminescence measurements were determined at the inoculation site. Twenty hours after inoculation of  $1 \times 10^6$  parasites, the bioluminescent signal was clearly detected at the inoculation site (Fig. 7A) and reached approximately  $1.8 \times 10^7$  Ph/s (Fig. 8A). Successful infection, defined as the development of lesions in ear dermis, was observed in 100% of the inoculated mice. In the groups of mice that were inoculated with  $1 \times 10^5$  *Leishmania* the infection rate was also 100% (not shown) but the bioluminescence was 15 times lower  $(1.2 \times 10^6 \text{ Ph/s}; \text{ Fig. 7B})$  than that in the group infected with the highest dose. In the last group of mice that were injected with  $1 \times 10^4$  parasites the bioluminescence was not detectable and not significantly different of the background (4 × 105 Ph/s; not shown). Eight days after inoculation of 1 × 10<sup>6</sup> parasites, the bioluminescent signal remained stable (Fig. 8) and reached  $2 \times 10^8$  Ph/s at day 16 after inoculation (Fig. 8A).

### **Discussion**

Deciphering biological processes in the context of tissues, systems requires the development of non-invasive imaging modalities such as *in vivo* bioluminescence imaging





**Fig. 8.** Monitoring the level of bioluminescence activity on ear after intradermal inoculation of 10<sup>6</sup> luciferase-transfected *Leishmania amazonensis*. Images were acquired with CCD camera over a 16 day time course and (A) quantification of luminescence in the ear dermis at day 1, 8 and 16 was monitored by detecting total photon emission per second. The background obtained for the same area was subtracted to the measured values. (B) One image of infected mouse at day 8 is displayed as pseudo-colour image, with variation in colour representing light intensities as described in previous figure. Colour scale next to image codes for signal intensity. This image is superimposed on the reference image for localization. On right panel, the three dimensional distribution and relative amplitude of luminescence are shown at the inoculation point. The highest line, corresponding to the red area, represents the most intense light emission.

(Greer and Szalay, 2002; Doyle *et al.*, 2004). In the present study we aimed to detect and track bioluminescent parasitic microorganisms through the course of the parasite/host interactions using the firefly luciferase as the reporter gene as well as its substrate the luciferin. The stable integration of the luciferase reporter gene into the ribosomal locus of *L. amazonensis* represents a valuable tool for (i) parasite targeted-drug discovery and testing, and (ii) assessing the discrete phases occurring in the different hosts they rely on.

Setting-up of standard curves for each environmental condition: the example of parasites displaying different developmental stages

Leishmania spp. display different developmental stages

that assess their reactivity to the host signals they rely on: while they develop as extracellular flagellated promastigotes in the insect vector, they develop as an obligate intracellular amastigotes that survive and multiply within the mammalian host. The stable integration of the luciferase reporter gene into the ribosomal locus of L. amazonensis represents a valuable tool for assessing phenotypes related to parasite features in laboratory mice when long periods of growth without selection is required. The measure of luciferase activity was compared and the level of sensitivity determined using recombinant parasites at different stages of differentiation and purified from culture or mouse lesions. It was possible to detect 10 times more luciferase-expressing amastigotes parasites extracted from cultured bone marrow-derived mouse macrophages, compared to amastigotes purified from mouse cutaneous lesions. Similarly when luciferase-expressing amastigotes derived from cultured macrophages, are differentiating into promastigotes, for the same number of parasites, we observed a 30-fold increase in luciferase activity in the latter. Several factors can contribute to lower luciferase activity in amastigotes and intracellular parasites such as their slower metabolism, different environmental conditions such as medium, protease inhibitors. Whatever the reasons are, the correlation between luciferase activity and parasite number was excellent both in culture and in infected macrophages or mouse tissues: as a result, standard curves for each condition have been set up and the level of sensitivity is relevant for most purposes and reproducible within one defined condition. The minimum number of parasites (amastigote stage) detected by their bioluminescent phenotypes was 1000 and in this study we show that the linear relationship of bioluminescence versus number of parasites was shown to give a dynamic measurement range of over four logs. Furthermore, it is possible to use luciferase activity derived from a standard curve established with purified amastigotes to determine parasite burdens in macrophages or in mouse tissues.

Setting-up of relevant readout assays for parasite targeted-drug discovery and testing

For a number of intracellular pathogenic microorganisms, the use of reporter genes has facilitated antimicrobial drug discovery and testing (Jacobs *et al.*, 1993; Buckner *et al.*, 1996; Collins *et al.*, 1998; Deb *et al.*, 2000; Chan *et al.*, 2003). Recently, models of axenic *Leishmania* amastigotes have been introduced and proposed for drug testing, instead of promastigotes (Callahan *et al.*, 1997; Ephros *et al.*, 1999; Sereno *et al.*, 2001). Still these axenic amastigotes differ from the amastigotes that either develop in unactivated/de-activated macrophages (Callahan *et al.*, 1997; Sereno *et al.*, 1998; Sereno *et al.*, 2001) or that persist in asymptomatic cutaneous tissue.

© 2004 Blackwell Publishing Ltd, Cellular Microbiology, 7, 383-392

In this study the activities of anti-leishmanial drugs against Leishmania sequestered in mouse bone marrowderived macrophages were evaluated by direct assay of the inhibition of luminescence. In order to validate the system this latter measure was compared with their effect on the evaluation of amastigote number in macrophage. The major difference between these two methods is that luciferase method measures the total number of parasites present while the staining method is based on the regression of parasitophorous vacuole (Fakhfakh et al., 2003) and the overall decrease in parasite number. Inhibition of luminescence correlated well with the leishmanicidal activity of various drugs tested. This level of this inhibition at the same concentration of drug was always similar to the inhibition of parasitophorous vacuole. In our conditions the level of sensitivity was sufficient to detect the presence of home parasite per 10 macrophages in a 96-well plate containing 4 × 10<sup>4</sup> host cells. Thus, luciferase-expressing parasites offer many advantages over alternative drug screening methods. This experimental system has merits, as it is relatively fast and inexpensive, and has potential for automation and high throughput drug screening. It will allow the rapid identification of candidate compounds that are suitable for further testing on amastigotes. In conclusion, from the data reported in this paper, it seems likely that recombinant Leishmania expressing reporter gene might be reliably and confidently used for the screening of new compounds active at least on DNA-replicating amastigotes that develop in mouse macrophages.

Imaging of bioluminescent parasites in mice: one way to visualize features of parasitism?

The next challenge was to dispose a direct assessment of the features of parasitism in real time. Does this technique provide both spatial and sequential information about the features of parasitism within the tissue context? Because of its non-destructive and non-invasive nature the imaging procedure can be performed repeatedly, allowing each animal to be used as its own control over time, overcoming the problem of animal-animal variations. Moreover, these in vivo methods do not require the removal of tissue and consequently the loss of contextual influences of the living animals (Contag et al., 1995; Francis et al., 2000; 2001; Rocchetta et al., 2001; Luker et al., 2002). More and more studies (Contag et al., 1995; Francis et al., 2000; 2001; Rocchetta et al., 2001; Luker et al., 2002) indicate not only that in vivo bioluminescent imaging is feasible but also that it can be quantitative and providing more information in less time than conventional assays. However, as far as the eukaryotic parasites are concerned, detection devices to visualize and quantify optical signal in living cells are still limited to in vitro or ex vivo analyses. Thus, Leishmania, transfected with reporter

genes, has been used in cell culture (Misslitz et al., 2000; Roy et al., 2000) and in ex vivo analyses after expression in animals (Misslitz et al., 2000; Roy et al., 2000; Chan et al., 2003; Sorensen et al., 2003). The present study is the first demonstration of detection and monitoring light emission in a murine model of parasitical infection. First, it was established that (i) the features of the parasites that have been transfected are not different from the nonluciferase expressing ones on the basis of the onset and the development of the lesion, and (ii) the intensity of the bioluminescence measured from the living mice correlated well with the parasite burden subsequently determined by standard protocols. Second, in the present study we observe that photon count imaging and the estimate of the cultivatable parasites increase in the same manner. Consequently, the bioluminescence could be used to monitor the parasitic load at the site of inoculation. In addition, following the intradermal delivery of low doses of invasive metacyclic promastigotes (Courret et al., 2003), the onset of the lesion in BALB/c mice was shown to assess the presence of doses as low as 10<sup>4</sup> parasites. The non-invasive approach we describe in the present paper could be further coupled to the use of mouse hosts expressing relevant bioluminescent probes especially if the sensitivity of the equipment is improved (Doyle et al., 2004).

In conclusion, this study presents a new model for visualizing the different discrete phases that Leishmania establishes in one of its hosts, the laboratory mice. The method described here is rapid, uses fewer animals and enables in vivo monitoring of the parasite burden. Additionally, this model may be useful for the study of pathogenesis and healing processes in mice, allowing to dissect one essential feature of parasitism, namely the parasite persistence, and its transmissible properties to the insect vector.

## **Experimental procedures**

Mice

We obtained 2 to 4 month old female BALB/c from Charles River (St Germain-sur-l'Arbresle, France). All the animal were handled under institutional guidelines of the Central Animal Facility at Institut Pasteur (Paris, France; http://www.pasteur.fr/recherche/ unites/animalerie/guidelin.html).

## Generation of bioluminescent Leishmania

A 1.66 kbp of firefly luciferase coding region was cut by Ncol/ Eagl from pGL3 basic (Promega, Madison, WI) and subsequently cloned into Leishmania expression vector pF4× 1.HYG (Jenabioscience, Jena, Germany) with marker gene for selection with Hygromycin B (Cayla, Toulouse, France) and previously cut by Ncol/Notl. In this construct 18 s rRNA flanked the luciferase and HYG genes. Following linearization with Swal, luciferase and HYG was integrated into 18 s rRNA locus of the nuclear DNA of *Leishmania*.

Transfection were realized by electroporation with following conditions:  $25~\mu F$ , 1500~V, in 4 mm cuvette;  $3.75~kV~cm^{-1}$  (Goyard *et al.*, 2003). Following electroporation, cells were incubated 24 h in media without drug, and plated on semisolid media containing  $100~\mu g~ml^{-1}$  of hygromycin B.

## Parasite preparation and inoculation in mice

Wild-type or transfected-L. amazonensis strain LV79 (MPRO/BR/ 1972/M1841) amastigotes were prepared from BALB/c mice (Antoine et al., 1989). Briefly, the promastigote form was grown at 26°C in M199 media supplemented with 10% FBS, 25 mM Hepes pH 6.9, 12 mM NaHCO<sub>3</sub>, 1 mM glutamine, 1× RPMI 1640 vitamin mix, 10 μM folic acid, 100 μM adenosine, 7.6 mM haemin. 50 U ml<sup>-1</sup> penicillin, and 50 ug ml<sup>-1</sup> streptomycin (Goyard et al., 2003). Infective-stage metacyclic promastigotes were isolated from stationary phase cultures (6 days old) by gradient of Ficoll as previously described (Spath and Beverley. 2001). Various doses (10<sup>4</sup>–10<sup>6</sup>) of metacyclic promastigotes per 10 µl of Dulbecco's PBS were injected in the ear dermis of BALB/ c mice. Lesion size was measured using a direct reading Vernier caliper (Thomas) and expressed as the difference in thickness between the inoculated ear and the non-inoculated contra-lateral ear.

#### Quantification of parasites

We estimated the number of parasites present in parasite-loaded ears and the draining lymph nodes, as previously described (Leclercq et al., 1996). The two sheets of the infected ears were separated, cut into small pieces and ground in HOSMEM-II culture medium, using a glass tissue homogenizer. Lymph nodes were removed, mechanically dissociated and the number of cells they contained was determined. The tissue/organ homogenates were serially diluted in HOSMEM-II culture medium and then dispensed into 96-well plates containing semisolid agar (Bacto-Agar, Difco, Detroit, MI) supplemented with 10% sterile heparintreated rabbit blood. The plates were incubated for 10 days and then each well was examined and classified as positive or negative according to whether or not viable promastigotes were present. Limiting dilution analysis was then applied to the data to estimate the number of viable parasites, expressed in LDAU. Statistical analysis of the results was based on the maximal likelihood method (Taswell, 1981; Strijbosch et al., 1987).

# Macrophage infections and parasites drug susceptibility assay

Macrophages were obtained from BALB/c mice by *in vitro* differentiation of bone marrow precursor cells in bacteriologic Petri dishes (Greiner, Nurtingen, FRG). Cells were cultured in RPMI 1640 medium (Seromed, Berlin, Germany) supplemented with 20% (v/v) heat-inactivated fetal calf serum (Dutscher, Brumath, France), 50 U ml $^{-1}$  penicillin, 50  $\mu g$  ml $^{-1}$  streptomycin and 20% L-929 fibroblast-conditioned medium. After 5 days at 37°C in a 5% CO $_2$ /95% air atmosphere, cells were washed in PBS and removed by treatment for 20 min with Ca $^{2+}$ , Mg $^{2+}$ -free PBS con-

taining 2 mg ml<sup>-1</sup> glucose (Prina et al., 1993). Recovered macrophages were further incubated in culture medium containing only 3% of conditioned medium either in 24- or 96-well plates or 35 mm culture dishes. Twenty-four hours after re-plating, macrophages were infected at a stationary phase promastigote-host cell ratio of 8:1 and then incubated at 34°C for 2 h before being washed with Dulbecco's phosphate-buffered saline (PBS, Seromed) to remove free parasites. Following 1 day incubation to allow differentiation in amastigotes, drug was added to the appropriate wells diluted in RPMI with 20% FBS and cells were incubated at 34°C for a further incubation of 24 h. For guinolines, compounds 22, 42 were used and the stock solutions were prepared in DMSO as described previously (Fakhfakh et al., 2003). Macrophages were either fixed and stained with Giemsa to assess the level of infection by light microscopy or directly treated for the quantification of bioluminescence. The percentage of amastigotes-loaded macrophages and the number of parasites per macrophages were determined by looking at least 200 macrophages in two independent wells.

# Extraction of intracellular parasites from macrophages and their further processing

Macrophages on 35 mm tissue cultures dishes were washed with PBS and then lysed in CHAPS lysis buffer (20 mM Tris-HCl buffer, pH 7.4, containing 130 mM NaCl and 0.3% (w/v) CHAPS (3[(3-chloroamidopropyl)dimethylammonio]-1-propane sulphonate, Sigma) (Courret *et al.*, 2001).

After 40 min at 4°C, macrophages were scraped off the dish using a rubber policeman. Recovered cells were collected and intracellular parasites were purified through a 45%/90% Percoll gradient (Pharmacia, Uppsala, Sweden). Viable parasites were washed and counted and prepared for luciferase assay.

# Estimation of luciferase content in transfected-Leishmania promastigotes and in infected macrophages

Parasites were washed in PBS and transferred in 96-well white culture plate (Dynex Technologies, Chantilly, France) to prevent cross-talk between wells during detection. Various doses of parasites were incubated with the luciferase assay system kit (Roche Molecular Biochemicals; Mannhein, Germany) for 5 min at room temperature. During this time complete cells lysis occurred and the light reaction started. Then plates were transferred to the Wallac MicroBeta Trilux scintillation counter (PerkinElmer Life Sciences, Boston, MA) used as luminometer and the light reaction was measured for 1 s. The luminescence units are luminescence counts per second (LCPS). In measuring luminescence the lower of the photomultiplier tubes (PMTs) is turned off and measurement is accomplished by the upper PMT only. Untranfected parasites were used as negative control. Leishmania-infected macrophages were washed with PBS. For measurement of bioluminescence, 100 µl of PBS was added to each well of macrophages and mixed with 100 µl of luciferase reagent.

# In vivo bioluminescence imaging of luciferase-expressing Leishmania

The mice were anesthetized by intraperitoneal administration of

ketamine (120 mg kg-1; Imalgène 1000, Mérial, France) and xylazine (4 mg kg<sup>-1</sup>; Rompun 2%, Bayer, Leverkusen, Germany). A solution of the luciferase substrate luciferin (150 mg kg<sup>-1</sup>; D-Luciferin potassium salt, Xenogen, California) was mixed with the anaesthetics just before use. A greyscale reference image of animals was generated and mice were imaged for a maximum of 5 min at a number of time points post-infection using a NightOwl cabinet with charge couple device (CCD) camera (Berthold Technologies, Bad Wildbad, Germany). Total photon emission from selected and defined areas within the images of each mouse was quantified with the WinLight32 software (Berthold). The photon signal from the ear was presented as a pseudocolour image representing light intensity (red most and blue least intense). This image was superimposed on the reference image for orientation.

#### **Acknowledgements**

The authors thank Frédérique Gautier for assistance with the Night Owl (Berthold Technologies) and Spencer Shorte for the warm welcome in the Plate-Forme d'Imagerie Dynamique (PFID) at the Institut Pasteur. The authors are extremely grateful to Alain Fournet for kindly providing quinolines. We also thank Eric Prina and Hélène Saklani for support and valuable advice. This work was supported by the Institut Pasteur and appel d'offres CNRS 'Puces à ADN' to Geneviève Milon together with M. Pagès/P. Glaser.

#### References

- Antoine, J.-C., Jouanne, C., and Ryter, A. (1989) Megasomes as the targets of leucine methyl ester in Leishmania amazonensis amastigotes. Parasitology 99: 1-9.
- Antoine, J.-C., Jouanne, C., Lang, T., Prina, E., de Chastellier, C., and Frehel, C. (1991) Localization of major histocompatibility complex class II molecules in phagolysosomes of murine macrophages infected with Leishmania amazonensis. Infect Immun 59: 764-775.
- Berman, J.D. (1984) Leishmania tropica: quantitation of in vitro activity of antileishmanial agents by Giemsa staining, viability, and 3H-formycin B incorporation. J Parasitol 70: 561-562.
- Blackwell, J.M. (1996) Genetic susceptibility to leishmanial infections: studies in mice and man. Parasitology 112: S67-S74.
- Buckner, F.S., Verlinde, C.L., La Flamme, A.C., and Van Voorhis, W.C. (1996) Efficient technique for screening drugs for activity against Trypanosoma cruzi using parasites expressing beta-galactosidase. Antimicrob Agents Chemother 40: 2592-2597.
- Buffet, P.A., Sulahian, A., Garin, Y.J., Nassar, N., and Derouin, F. (1995) Culture microtitration: a sensitive method for quantifying Leishmania infantum in tissues of infected mice. Antimicrob Agents Chemother 39: 2167-2168.
- Callahan, H.L., Portal, A.C., Devereaux, R., and Grogl, M. (1997) An axenic amastigote system for drug screening. Antimicrob Agents Chemother 41: 818-822.
- Chan, M.M.Y., Bulinski, J.C., Chang, K.P., and Fong, D. (2003) A microplate assay for Leishmania amazonensis promastigotes expressing multimeric green fluorescent protein. Parasitol Res 89: 266-271.
- Collins, L.A., Torrero, M.N., and Franzblau, S.G. (1998)

- Green fluorescent protein reporter microplate assay for high-throughput screening of compounds against Mycobacterium tuberculosis. Antimicrob Agents Chemother 42: 344-347.
- Contag, C.H., Contag, P.R., Mullins, J.I., Spilman, S.D., Stevenson, D.K., and Benaron, D.A. (1995) Photonic detection of bacterial pathogens in living hosts. Mol Microbiol 18: 593-603.
- Courret, N., Fréhel, C., Prina, E., Lang, T., and Antoine, J.-C. (2001) Kinetics of the intracellular differentiation of Leishmania amazonensis and internalization of host MHC molecules by the intermediate parasite stages. Parasitology 122: 263-279.
- Courret, N., Lang, T., Milon, G., and Antoine, J.C. (2003) Intradermal inoculations of low doses of Leishmania major and Leishmania amazonensis metacyclic promastigotes induce different immunoparasitic processes and status of protection in BALB/c mice. Int J Parasitol 33: 1373-1383.
- Deb, D.K., Srivastava, K.K., Srivastava, R., and Srivastava, B.S. (2000) Bioluminescent Mycobacterium aurum expressing firefly luciferase for rapid and high throughput screening of antimycobacterial drugs in vitro and in infected macrophages. Biochem Bioph Res Co 279: 457-461.
- Doyle, T.C., Burns, S.M., and Contag, C.H. (2004) In vivo bioluminescence imaging for integrated studies of infection. Cellular Microbiol 6: 303-317.
- Edinger, M., Hoffmann, P., Contag, C.H., and Negrin, R.S. (2003) Evaluation of effector cell fate and function by in vivo bioluminescence imaging. Methods 31: 172-179.
- Ephros, M., Bitnun, A., Shaked, P., Waldman, E., and Zilberstein, D. (1999) Stage-specific activity of pentavalent antimony against Leishmania donovani axenic amastigotes. Antimicrob Agents Chemother 43: 278-282.
- Fakhfakh, M.A., Fournet, A., Prina, E., Mouscadet, J.F., Franck, X., Hocquemiller, R., and Figadere, B. (2003) Synthesis and biological evaluation of substituted quinolines: Potential treatment of protozoal and retroviral co-infections. Bioora Med Chem 11: 5013-5023.
- Francis, K.P., Joh, D., Bellinger-Kawahara, C., Hawkinson, M.J., Purchio, T.F., and Contag, P.R. (2000) Monitoring bioluminescent Staphylococcus aureus infections in living mice using a novel luxABCDE construct. Infect Immun 68: 3594-3600.
- Francis, K.P., Yu, J., Bellinger-Kawahara, C., Joh, D., Hawkinson, M.J., Xiao, G., et al. (2001) Visualizing pneumococcal infections in the lungs of live mice using bioluminescent Streptococcus pneumoniae transformed with a novel grampositive lux transposon. Infect Immun 69: 3350-3358.
- Goyard, S., Segawa, H., Gordon, J., Showalter, M., Duncan, R., Turco, S.J., and Beverley, S.M. (2003) An in vitro system for developmental and genetic studies of \*Leishmania\* donovani phosphoglycans. Mol Biochem Parasitol 130:
- Greer, L.F., and Szalay, A.A. (2002) Imaging of light emission from the expression of luciferases in living cells and organisms: a review. Luminescence 17: 43-74.
- Handman, E. (2001) Leishmaniasis: Current Status of Vaccine Development. Clin Microbiol Rev 14: 229-243.
- Jacobs, W.R., Barletta, R.G., Udani, R., Chan, J., Kalkut, G., Sosne, G., et al. (1993) Rapid assessment of drug susceptibilities of mycobacterium-tuberculosis by means of Luciferase Reporter Phages. Science 260: 819-822.
- Kadurugamuwa, J.L., Sin, L., Albert, E., Yu, J., Francis, K., DeBoer, M., et al. (2003) Direct continuous method for

- monitoring biofilm infection in a mouse model. Infect Immun 71: 882-890.
- Katakura, K., Kawazu, S., Naya, T., Nagakura, K., Ito, M., Aikawa, M., et al. (1998) Diagnosis of kala-azar by nested PCR based on amplification of the Leishmania mini-exon gene. J Clin Microbiol 36: 2173-2177.
- Launois, P., Louis, J.A., and Milon, G. (1997) The fate and persistence of Leishmania major in mice of different genetic backgrounds: an example of exploitation of the immune system by intracellular parasites. Parasitology 115: S25-S32.
- Leclercq, V., Lebastard, M., Belkaid, Y., Louis, J., and Milon, G. (1996) The outcome of the parasitic process initiated by Leishmania infantum in laboratory mice: a tissuedependent pattern controlled by the Lsh and MHC loci. J Immunol 157: 4537-4545.
- Loimaranta, V., Tenovuo, J., Koivisto, L., and Karp, M. (1998) Generation of bioluminescent Streptococcus mutans and its usage in rapid analysis of the efficacy of antimicrobial compounds. Antimicrob Agents Chemother 42: 1906-1910.
- Luker, G.D., Bardill, J.P., Prior, J.L., Pica, C.M., Piwnica-Worms, D., and Leib, D.A. (2002) Noninvasive bioluminescence imaging of herpes simplex virus type 1 infection and therapy in living mice. J Virol 76: 12149-12161.
- Lugue-Ortega, J.R., Rivero-Lezcano, O.M., Croft, S.L., and Rivas, L. (2001) In vivo monitoring of intracellular ATP levels in Leishmania donovani promastigotes as a rapid method to screen drugs targeting bioenergetic metabolism. Antimicrob Agents Chemother 45: 1121-1125.
- Misslitz, A., Mottram, J.C., Overath, P., and Aebischer, T. (2000) Targeted integration into a rRNA locus results in uniform and high level expression of transgenes in Leishmania amastigotes. Mol Biochem Parasitol 107: 251-261.
- Nicolas, L., Milon, G., and Prina, E. (2002) Rapid differentiation of Old World Leishmania species by LightCycler polymerase chain reaction and melting curve analysis. J Microbiol Methods 51: 295-299.
- Noyes, H.A., Reyburn, H., Bailey, J.W., and Smith, D. (1998) A nested-PCR-based schizodeme method for identifying Leishmania kinetoplast minicircle classes directly from clinical samples and its application to the study of the epidemiology of Leishmania tropica in Pakistan. [see comments] J Clin Microbiol 36: 2877-2881.

- Prina, E., Jouanne, C., de Souza Lao, S., Szabo, A., Guillet, J.G., and Antoine, J.-C. (1993) Antigen presentation capacity of murine macrophages infected with Leishmania amazonensis amastigotes. J Immunol 151: 2050-
- Rocchetta, H.L., Boylan, C.J., Foley, J.W., Iversen, P.W., Letourneau, D.L., McMillian, C.L., et al. (2001) Validation of a noninvasive, real-time imaging technology using bioluminescent Escherichia coli in the neutropenic mouse thigh model of infection. Antimicrob Agents Chemother 45: 129-137.
- Roy, G., Dumas, C., Sereno, D., Wu, Y., Singh, A.K., Tremblay, M.J., et al. (2000) Episomal and stable expression of the luciferase reporter gene for quantifying Leishmania spp. infections in macrophages and in animal models. Mol Biochem Parasitol 110: 195-206.
- Sereno, D., Cavaleyra, M., Zemzoumi, K., Maguaire, S., Ouaissi, A., and Lemesre, J.L. (1998) Axenically grown amastigotes of Leishmania infantum used as an in vitro model to investigate the pentavalent antimony mode of action. Antimicrob Agents Chemother 42: 3097-3102.
- Sereno, D., Roy, G., Lemesre, J.L., Papadopoulou, B., and Ouellette, M. (2001) DNA transformation of Leishmania infantum axenic amastigotes and their use in drug screening. Antimicrob Agents Chemother 45: 1168-1173.
- Sorensen, M., Lippuner, C., Kaiser, T., Misslitz, A., Aebischer, T., and Bumann, D. (2003) Rapidly maturing red fluorescent protein variants with strongly enhanced brightness in bacteria. FEBS Lett 552: 110-114.
- Spath, G.F., and Beverley, S.M. (2001) A lipophosphoglycanindependent method for isolation of infective Leishmania metacyclic promastigotes by density gradient centrifugation. Exp Parasitol 99: 97-103.
- Strijbosch, L.W., Buurman, W.A., Does, R.J., Zinken, P.H., and Groenewegen, G. (1987) Limiting dilution assays. Experimental design and statistical analysis. J Immunol Methods 97: 133-140.
- Taswell, C. (1981) Limiting dilution assays for the determination of immunocompetent cell frequencies. I. Data analysis. J Immunol 126: 1614-1619.
- Titus, R.G., Marchand, M., Boon, T., and Louis, J.A. (1985) A limiting dilution assay for quantifying Leishmania major in tissues of infected mice. Parasite Immunol 7: 545-