

1 Indotecan (LMP400) and AM13-55: Two novel indenoisoquinolines show potential
2 for treating visceral leishmaniasis

3

4 Rafael Balaña-Fouce¹, Christopher F. Prada¹, José María Requena², Mark Cushman³,
 5 Yves Pommier⁴, Raquel Álvarez-Velilla¹, José Miguel Escudero-Martínez¹, Estefania
 6 Calvo-Álvarez¹, Yolanda Pérez-Pertejo¹, Rosa M. Reguera¹

7

8 ¹Departamento de Ciencias Biomédicas, Universidad de León, Campus de Vegazana
 9 s/n; 24071 León (SPAIN), Phone 34 987 291257; Telefax 34 987 291252

10 ²Departamento de Bioquímica y Biología Molecular, Centro de Biología Molecular
 11 “Severo Ochoa” c/ Nicolás Cabrera, 1, 28049 Madrid (SPAIN)

12 ³Department of Medicinal Chemistry and Molecular Pharmacology, College of
 13 Pharmacy, and the Purdue Center for Cancer Research, Purdue University, West
 14 Lafayette, IN 47907, USA

15 ⁴Laboratory of Molecular Pharmacology, Center for Cancer Research, National Cancer
 16 Institute, Building 37, Room 5068, 37 Convent Drive, Bethesda, Maryland 20892,
 17 (USA)

18 **Correspondence to:** Dr. Rafael Balaña-Fouce. Departamento de Ciencias Biomédicas,
 19 Universidad de León, Campus de Vegazana s/n; 24071 León (SPAIN), Phone 34 987
 20 291257; telefax 34 987 291252; e-mail: rbalf@unileon.es

21 **Running title:** Leishmanicide effect of indenoisoquinolines

22

23 **Keywords:** DNA topoisomerase IB, indenoisoquinolines; infrared protein; splenic
 24 explants; *Leishmania*

25

26 **Abbreviations:** NTD: Neglected Tropical Diseases; Top: DNA-topoisomerases;
 27 LiTopIB: Type IB *Leishmania infantum* Top; hTopIB: Type IB human Top; DALY:
 28 Disability-Adjusted Life Year; CPT: camptothecin; IFP1.4: Infrared Fluorescent Protein
 29 1.4; FCS: foetal calf serum

30 **ABSTRACT**

31 Visceral leishmaniasis is an emerging neglected tropical disease (NTD) caused by the
32 protozoan *Leishmania infantum* in the countries bordering the Mediterranean Basin.
33 Currently there is no effective vaccine against this disease, and the therapeutic approach
34 is based in toxic derivatives of Sb^V. Therefore, the discovery of new therapeutic targets
35 and the development of drugs designed to inhibit them is an extremely important
36 approach to fight this disease. DNA topoisomerases (Top) have been identified as
37 promising targets against leishmaniasis. These enzymes are involved in solving
38 topological problems generated during replication, transcription and recombination of
39 DNA. Unlike the mammalian host, type IB DNA topoisomerase (TopIB) from
40 *Leishmania* spp. is a unique bisubunit protein that makes it very interesting as a
41 selective drug target. In the present investigation, we studied the effect of two TopIB
42 poisons with indenoisoquinoline structure; Indotecan and AM13-55, on a murine
43 BALB/c model of infected splenocytes with *L. infantum*, comparing their effectiveness
44 with the clinically tested leishmanicide drug, paromomycin. Both compounds have
45 high selectivity indexes when compared with uninfected splenocytes. SDS/KCl-
46 precipitable DNA-protein complexes in *Leishmania* promastigotes and *in vitro* cleaving
47 assays confirmed that these drugs are Top poisons. The inhibitory potency of both
48 indenoisoquinolines on *L. infantum* recombinant TopIB was assessed *in vitro* showing
49 that indotecan was the most active compound preventing the relaxation of supercoiled
50 DNA. Experimental infections in susceptible BALB/c mice treated with 2.5 mg/kg body
51 weight/day once every other day for a total of fifteen days showed that indotecan
52 cleared more than 80% parasite burden of spleen and liver indicating promising activity
53 against visceral leishmaniasis.

54

55 **INTRODUCTION**

56 Neglected Tropical Diseases (NTD) produced by vector-borne protozoa are mostly
57 prevalent in underdeveloped and poor countries causing an estimated 4.1 million
58 disability-adjusted life years (DALYs) lost (16). However, in developed countries
59 where these diseases were unknown or had been eradicated for a long time, an
60 unexpected new scenario has appeared. Increased tourist exchange, soldiers deployed in
61 countries where these diseases are endemic, migratory fluxes and eventually climate
62 changes, along with other pandemics, are challenging the apparent safety of rich
63 populations (26,30). Since NTDs are prevalent in low-income countries, large
64 pharmaceutical companies have neglected R & D of new drugs. Therefore, old
65 compounds that are losing efficacy are still in use. These compounds have undesirable
66 toxic effects and their dosage schedules are complex and repetitive, which compromises
67 patient compliance (7).

68 Visceral leishmaniasis is a NTD increasingly affecting European countries by the
69 massive flows of immigrants from North Africa (18). Its etiologic agent *Leishmania*
70 *infantum* infects domestic dogs (canine leishmaniasis) that act as reservoir, and it is
71 transmitted to humans through sandfly bites (20). For decades, the first-line treatment of
72 this disease was the old-fashioned and toxic drugs derived from pentavalent antimony
73 (Sb^V) (1). Recently, safer and more effective drugs such as amphotericin B, miltefosine
74 and paromomycin are substituting Sb^V , despite the fact that they are not devoid of
75 undesirable side effects and cannot be administered during pregnancy (8,9). Based on
76 these reasons, the search for new compounds against this disease is very much needed.
77 Since the beginning of this century, DNA topoisomerase IB (TopIB) has been identified
78 as a potential target against *Leishmania* and other trypanosomatids (3,5,6). The choice
79 of this target is based on two main reasons: i) the enzyme has an increased expression
80 during the rapidly dividing cycle of the pathogen, in a similar way to tumour cells, and
81 especially ii) the pathogen's enzyme is structurally different from that of the host (35).

82 Our research group found that *L. donovani* TopIB (LdTopIB) was atypically composed
83 of two different subunits – each one encoded by different genes – that have to be
84 assembled in the pathogen to reconstitute the active enzyme. One of the subunits
85 contains the four amino acids of the active site, which is fully conserved from a
86 phylogenetic point of view (35). The other subunit contains the catalytic amino acid
87 (Tyr222), which acts by breaking one DNA strand by a specific nucleotide sequence

88 (13). All these features have also been described in the other protozoan-borne NTDs
89 pathogens: *Trypanosoma brucei* – the agent responsible for sleeping sickness in Africa
90 (11) – and *T. cruzi* – responsible for Chagas disease in South America (36). However,
91 despite the fact that these enzymes conserve their catalytic domains unchanged, they
92 display two non-conserved regions; one at the C-terminal end of the large protomer and
93 the other at the N-terminal end of the small protomer, which are extremely important in
94 sensitivity to topoisomerase poisons (12). These compounds act by stabilizing enzyme-
95 DNA complexes, preventing the religation step and eventually producing single-strand
96 breaks when they collide with the replication fork during DNA synthesis (22). The most
97 studied TopIB poison is camptothecin (CPT) and its derivatives, including topotecan
98 and the prodrug irinotecan, which are being used against certain tumours (32). Other
99 non-CPT TopIB poisons are indolocarbazoles, such as the DNA intercalating drug
100 rebeccamycin, and indenoisoquinolines, which were initially developed as antitumor
101 compounds with an improved ability to stabilize cleavage complexes (2,25).

102 There are several promising indenoisoquinolines studied against tumour processes and
103 more recently against experimental African trypanosomiasis (4). In this paper we show
104 for the first time the effect of two indenoisoquinolines, indotecan and AM13-55, on the
105 viability of free-living promastigotes and amastigote-infected murine splenocytes. For
106 this purpose, we used a modified strain of *L. infantum* transfected with the gene
107 encoding the infrared protein IFP1.4 from the extremophilic bacterium *Deinococcus*
108 *radiodurans* (31). This fluorescent pathogen strain retains its virulence unaffected and
109 allows studies of high performance in both free-living promastigotes and in *ex vivo*
110 spleen explants where parasite growth conditions resemble natural infection
111 environment. The results provided in this work demonstrate the high selectivity indexes
112 of both molecules *in vitro* and the promising therapeutic potential of indotecan against
113 visceral leishmaniasis.

114

115 **MATERIALS AND METHODS**

116 **Reagents and culture media**

117 *Pyrococcus furiosus* (*Pfu*), klenow polymerases and restriction enzymes were acquired
118 from Roche (Roche Farma SA, Spain) and GE Healthcare (Spain). T4 DNA ligase was
119 obtained from Stratagene (La Jolla, CA, USA). Cell culture media were purchased from

120 Sigma-Aldrich (Spain). Indenoisoquinolines indotecan and AM13-55 (Fig. 1) were
 121 kindly provided by Dr. Mark Cushman, Dept. of Medicinal Chemistry and Molecular
 122 Pharmacology, Purdue University (USA). Primers for PCR amplification were from
 123 Sigma Genosys (UK).

124

125 **Generation of an infrared fluorescent *L. infantum* strain**

126 *L. infantum* promastigotes (BCN150 strain) were obtained by Dr. J. M. Requena
 127 (Centro de Biología Molecular “Severo Ochoa”, Madrid Spain). The 987 bp IFP1.4
 128 coding region – an improved monomeric infrared fluorescent protein derived from
 129 *Deinococcus radiodurans* (31) – was amplified by PCR from IFP1.4_pcDNA3 vector –
 130 a gift from Dr. Roger Y. Tsien (Department of Pharmacology; Department of Chemistry
 131 and Biochemistry, University of California, San Diego, USA) – using the forward
 132 primer 5’CCGCTCGAGCCATGGCCACCATGGCTCGGGACCCTCTGCC3’ and the
 133 reverse primer 5’ATAAGAATGCGGCCGCTCATTTATACAGCTCGTCCATTCC3’.
 134 The amplified fragment was digested with *Bgl*II and *Not*I and assembled into the
 135 expression vector pLEXSY-2-HYG (Jena Bioscience GmbH; Germany), which was
 136 previously digested with the same enzymes. The resultant plasmid was named
 137 pLEXSY-IFP1.4. Parasites were electroporated with the large *Swa*I targeting fragment
 138 from pLEXSY-IFP1.4 and selected on semisolid media as previously described (27).
 139 Many clonal lines were obtained, and correct integration into the 18s rRNA locus was
 140 confirmed by PCR analysis. The transformant strain was routinely cultured at 26 °C in
 141 M199 medium supplemented with 25 mM HEPES pH 6.9, 10 mM glutamine, 7.6 mM
 142 hemin, 0.1 mM adenosine, 0.01 mM folic acid, 1x RPMI 1640 vitamin mix (Sigma),
 143 10% (v/v) heat-inactivated foetal calf serum (FCS) and antibiotic cocktail (50 U/ml
 144 penicillin, 50 µg/ml streptomycin).

145

146 ***Ex vivo* splenic explant cultures**

147 Balb/c mice were inoculated intraperitoneally with 10⁸ *L. infantum* purified metacyclic
 148 promastigotes. Briefly, infective promastigotes were isolated from stationary-phase
 149 culture by negative selection with peanut agglutinin (29). Five weeks post-infection
 150 spleens were aseptically dissected, washed in cold-PBS and placed in Petri dishes.
 151 Small pieces were obtained by using a scalpel. In order to obtain a single cell

suspension, tissue was incubated with 5 ml of 2 mg/ml collagenase D (Roche) prepared in buffer (10 mM Hepes pH 7.4, 150 mM NaCl, 5 mM KCl, 1 mM MgCl₂, 1.8 mM CaCl₂) for 20 min at 37 °C. Cell suspension containing traces of spleen mass was gently passed through a 100 µm cell strainer to remove tissue fragments (21). Splenocytes were washed twice with PBS by centrifugation (500 x g for 7 min at 4 °C) and re-suspended in RPMI medium supplemented with 10% FCS, 1 mM sodium pyruvate, 1x RPMI vitamins, 10 mM Hepes, and 50 U/ml penicillin and 50 µg/ml streptomycin at 37 °C and 5% CO₂. Cells were counted and diluted at different cell densities. Cells were seeded until confluence and different concentrations of the studied indenoisoquinolines were administered to the explants for 48 h. The viability of infecting amastigotes was assessed by registering the fluorescence emission at 708 nm in an Odyssey (Li-Cor) infrared imaging system. The cytotoxicity of the drugs was assessed on uninfected *ex vivo* explants after 48 h incubation, using the Alamar Blue staining method according to manufacturer's recommendations.

166

167 **Recombinant *Leishmania infantum* TopIB (LiTopIB)**

168 Cloning of *LiTopIB* ORFs (encoding large and small subunits), expression and
 169 purification of the enzyme were carried out as previously described for LdTopIB (35).
 170 Briefly, a *Saccharomyces cerevisiae* EKY3 strain deficient in TopIB activity [*MAT α*
 171 *ura3-52 his3Δ200 leu2 Δ1 trp1 Δ63 top1 Δ:TRP1*], was transformed by the lithium
 172 acetate method (15) with the bicistronic pESC-URA vector, which carries both *LiTopIB*
 173 ORFs. Single colonies were incubated overnight in SC-uracil media with 2% dextrose
 174 (w/v). Since dextrose traces may interfere with the expression of the protein, yeast
 175 cultures were incubated in SC-uracil media supplemented with 2% raffinose (w/v) for
 176 24 h prior to a 6-h induction with 2% galactose (w/v). Yeasts were harvested, washed
 177 with cold 1 x TEEG buffer (50 mM Tris-HCl pH 7.4, 1 mM EDTA, 1 mM EGTA, 10%
 178 glycerol) and resuspended in 15 ml of the same buffer supplemented with 0.2 M KCl
 179 and a protease inhibitors cocktail (Roche Farma SA, Spain). The cells were subjected to
 180 one freeze/thaw cycle at 80 °C and lysed by vortexing in a glass bead-beater at 4 °C.
 181 Protein extracts were obtained by centrifugation at 15,000 x g for 45 min at 4 °C. Yeast
 182 extracts were sequentially precipitated with two increasing concentrations of
 183 ammonium sulfate (35 and 75%, respectively). The second precipitate supernatant was
 184 loaded onto a P-11 phosphocellulose column (Whatman International Ltd. England),

185 which was equilibrated following manufacturer's indications. LiTopIB protein was
186 eluted at 4 °C with a discontinuous gradient of KCl (0.2, 0.4, 0.6, 0.8 and 1 M) in TEEG
187 buffer supplemented with 0.1 mg/ml sodium bisulphite, 0.8 mg/ml sodium bisulfite and
188 the protease inhibitors cocktail. Active fractions were loaded onto a phenyl-sepharose
189 column (Sigma-Aldrich) with a discontinuous inverse gradient of ammonium sulfate (1,
190 0.8, 0.6, 0.4 and 0.2 M). Elution was performed with subsequent Centricon (Millipore®)
191 concentration.

192

193 **Supercoiled plasmid DNA TopIB-mediated cleavage assay**

194 The sensitivity of LiTopIB to indotecan, AM13-55 and CPT (used as control drug) was
195 assayed by DNA cleavage assays. pBluescript SK(-) phagemid DNA (pSK) was used as
196 nickable substrate. At least 100 units of purified LiTopIB were incubated with 0.5 µg of
197 pSK DNA in 10 mM Tris-HCl buffer pH 7.5, 5 mM MgCl₂, 5 mM DTT, 0.1 mM
198 EDTA, 15 µg/ml bovine serum albumin and 50 mM KCl. Different drug concentrations
199 in 1% (v/v) DMSO (20 µl reaction volume) were also added. Following incubation for 4
200 min at 25 °C, reactions were then stopped with up to 1% (w/v) SDS and incubated in the
201 presence of 1 mg/ml proteinase K for one extra hour at 37 °C. Subsequently samples
202 were extracted with one volume of phenol-chloroform and loaded onto a 1% agarose gel
203 containing ethidium bromide to a final concentration of 40 pg/µl (17). The gel was run
204 for 16 h at 4 V/cm and the images of cleavage products were acquired with a G-BOX
205 (Syngene UK)

206

207 **DNA relaxation assays**

208 TopIB activity was assayed by the relaxation of negatively supercoiled plasmid DNA.
209 One unit of recombinant LiTopIB was incubated with the corresponding drug for 15
210 min at 4 °C. Then, the reaction mixture containing 0.5 µg of supercoiled pSK DNA, 10
211 mM Tris-HCl buffer pH 7.5, 5 mM MgCl₂, 0.1 mM EDTA, 15 µg/mL bovine serum
212 albumin, 50 mM KCl, in a total volume of 20 µl was added. Reaction mixtures were
213 incubated for 4 min at 25 °C. Enzyme reactions were stopped by the addition of up to
214 1% SDS (w/v) (final concentration) and digested with 1 mg/ml proteinase K at 37 °C
215 during one extra hour to remove the protein that remained linked to DNA fragments.
216 The extent of plasmid DNA relaxation was assessed in 1% agarose gels by

217 electrophoresis in 0.1 M Tris borate EDTA buffer (pH 8.0) at 2 V/cm for 16 h. Gels
218 were visualized with UV illumination after ethidium bromide (0.5 µg/ml) staining. A
219 further electrophoresis was run in the presence of 0.1 µg/ml ethidium bromide, in order
220 to separate nicked DNA from relaxed topoisomers (33).

221

222 **SDS-KCl precipitation assay**

223 For indenoisoquinoline-induced protein-DNA complex determination, *L. infantum*
224 promastigotes, previously labeled for 24 h with 0.5 µCi/ml [2-¹⁴C] thymidine, were
225 exposed to different concentrations of CPT and indenoisoquinolines for 30 min (3).
226 Cells were pelleted and lysed by incubation at 60 °C for 10 min in the presence of
227 1.25% (w/v) SDS; 0.4 mg/ml salmon sperm DNA and 5 mM EDTA. After the addition
228 of 65 mM KCl, the reaction mixture was incubated on ice for 60 min. The precipitate
229 was harvested by filtering through glass fiber paper (GF/C; Brandel Inc. MD, USA),
230 pre-wetted with wash buffer (10 mM Tris-HCl, pH 8, 1 mM EDTA, 100 mM KCl),
231 washed (three times with 4 ml of wash buffer), and air-dried. After spotting 50 µl of the
232 same density of labeled cell suspensions onto filter paper and precipitating nucleic acids
233 using 5% trichloroacetic acid (v/v), total incorporation of radioactivity into DNA was
234 measured by a liquid scintillation counter (Beckman LS600A). All assays included
235 vehicle drug as control. The formation of DNA fragments, as a percentage of total
236 labeled DNA, was calculated as follows: [(dpm in SDS-KCl drug – dpm in SDS-KCl
237 solvent)/(dpm total incorporation)] x 100.

238

239 **Thymidine incorporation assay**

240 *Leishmania infantum* promastigotes (10 ml; 5×10^6 cells/ml) were incubated with [2-
241 ¹⁴C] thymidine (0.5 µCi/ml) in the presence of drug or solvent in thymidine-free
242 complete M199 medium. Five hundred µl aliquots were removed after 2, 4, 8, 10 and 24
243 hours and loaded onto glass fiber paper (GF/C; Brandel Inc. MD, USA). ¹⁴C-labeled-
244 DNA was precipitated using ice-cold 5% (w/v) trichloroacetic acid and sequentially
245 washed with PBS, 90% (v/v) ethanol and 70% (v/v) ethanol. Radioactivity was
246 quantified by liquid scintillation. CPT was used as positive control.

247

248 **Animals**

249 Female BALB/c mice (age 4–6 weeks) were purchased from Harlan Ibérica
250 Laboratories (Spain). All animal procedures were approved by the University of Leon
251 subcommittee on Research Animal Care.

252 Infective-stage metacyclic *L. infantum* promastigotes were isolated from stationary
253 phase cultures (6-day old) by negative selection with peanut agglutinin (29). Mice were
254 infected with 10^7 metacyclic parasites intravenously through the tail vein. Fourteen days
255 later, once every two-days for fifteen days, mice were administered a solution of
256 indotecan or AM13-55 indenoisoquinolines diluted in DMSO (equivalent to a dose of
257 2.5 mg/kg of body weight per injection) intraperitoneally for a total of eight doses. Five
258 days after the last treatment, the mice were euthanized, organs were removed and
259 parasite loads were determined by limiting dilution assay (34). Briefly, a suspension of
260 20 mg of either spleen or liver was prepared by grinding the tissue in 1 ml of
261 Schneider's medium (Gibco BRL) containing 20% (v/v) heat-inactivated FCS in the
262 presence of streptomycin and penicillin. This suspension was further diluted to reach a
263 final concentration of 1 mg/ml. The parasite burden was determined by the limit dilution
264 method after a 10 day period of parasite growth. Parasite burden was expressed as the
265 number of free-living promastigotes per gram of wet tissue and compared to the non-
266 treated group.

267

268 **RESULTS AND DISCUSSION**

269 Recent studies showed that the CPT derivatives topotecan, SN38, and especially
270 gimatecan, are powerful leishmanicide agents. *In vitro* assays demonstrated that they act
271 by trapping both DNA and LiTopIB in reversible ternary complexes, producing
272 precipitable SDS/KCl material when free-living promastigotes were incubated in the
273 presence of these drugs (Prada et al., in press) (23). These compounds had good
274 selectivity indexes when their cytotoxicity on mammalian cells (un-infected
275 splenocytes) was compared to that from isolated *ex vivo* splenic explants of BALB/c
276 mice infected with *L. infantum*. However, none of these compounds were more selective
277 than miltefosine, the only effective oral treatment for visceral and cutaneous
278 leishmaniasis (9). The use of a model of mouse splenocytes infected with a *L. infantum*
279 strain expressing IFP1.4 permits high-throughput screening of compounds under

conditions that resemble those found in the animal, including the complete range of immune host cells, infected macrophages and fibroblasts (28).

Screening of indenoisoquinolinic compounds

The initial screening of new indenoisoquinolinic compounds was performed on free-living promastigotes; IRF1.4-*L. infantum* (BCN150) strain. To quantify cytotoxicity of the test compounds we utilized peritoneal BALB/c mouse macrophages. We first excluded compounds, that fell below an arbitrary cytotoxicity threshold in the peritoneal macrophage cultures ($IC_{50} < 10 \mu M$) (data not shown), and after exclusion of nine indenoisoquinolinic compounds (Fig. S1), only two new drugs were kept for further validation studies (Fig 1).

For these two compounds, anti-leishmanial activity (EC_{50}) was determined in the *ex vivo* splenic explant system. Comparison of these EC_{50} values with cytotoxicity (IC_{50}) values in the un-infected *ex vivo* splenic explant allowed determination of an *in vitro* selectivity index (SI: IC_{50}/EC_{50}) (Table 1). As positive control we have used both paromomycin and CPT. Paromomycin has demonstrated a good selectivity index in the same system (Prada *et al.*, in press (23)).

CPT and both indenoisoquinolines (indotecan and AM13-55) were much more effective in preventing the proliferation of promastigotes and development of infection in splenocytes than paromomycin, an aminoglycoside antibiotic used in clinical practice against human leishmaniasis. The IC_{50} of paromomycin was estimated to be $9.20 \mu M$ for mouse *ex vivo* splenic explants infected with *L. infantum*, whereas the IC_{50} value was $0.1 \mu M$ for indotecan and AM13-55. Only CPT was more effective ($IC_{50} = 0.03 \mu M$) at killing the parasites. To estimate the cytotoxicity on mammalian cells and therefore, to determine selectivity indexes of the studied compounds, we prepared non-infected *ex vivo* splenic explant. These cells were exposed to different concentration ranges of the drugs for 48 h and the viability was determined by the Alamar Blue method. Dose response curves were fitted by nonlinear analysis with the Sigma-Plot statistical package, showing that the least toxic compound was paromomycin ($IC_{50} = 15.70 mM$), followed by indotecan ($IC_{50} = 57.16 \mu M$) and AM13-55 ($IC_{50} = 48.37 \mu M$). Selectivity indexes of each compound were calculated as the ratio between the IC_{50} values on the uninfected explants system vs. the IC_{50} values on infected *ex vivo* splenic explants. Both

indenoisoquinolines have very high selectivity indexes; 571.6 for indotecan and 483.7 for AM13-55, which are very much higher than CPT-related compounds and miltefosine (23), but less than paromomycin (the safest compound tested with a selectivity index of 1706.5).

Indotecan and AM13-55 induce TopI-DNA covalent complexes

Figure 2A shows the induction of DNA cleavage complexes by indotecan and AM13-55 in the presence of TopI as tested in supercoiled plasmid DNA (pSK). Both indotecan and AM13-55 induce DNA cleavage complexes in a similar pattern to CPT, but with differences in their relative intensities. The indotecan cleavage complex showed the same strong intensity than CPT at 100 μ M, although the cleavage complex intensity is almost the same in the range of 0.1 to 10 μ M. On the other hand, the AM13-55 cleavage complex has lower intensities as CPT but the pattern resembles CPT, being proportional with drug concentration. We observed a dose-dependent increase in nicked DNA for the three compounds, thus confirming the TopIB poisoning nature of both indenoisoquinolines (2,4).

We also studied the potency to inhibit the DNA relaxation activity, comparing the human and leishmania TopIB enzymes. The LiTopIB enzyme was more sensitive to indotecan and AM13-55 than the human enzyme. Figure 2B (top panel) showed that Leishmania TopIB was already inhibited at 80 nM indotecan. However, 2.5 μ M were required to achieve the same effect on the human enzyme. AM13-55 (Fig 2B lower panel) was a little bit less efficient, inhibiting the *Leishmania* TopIB at 0.15 μ M, while the human enzyme was not inhibited until 2.5 μ M.

These results suggest that indenoisoquinolines indotecan and AM13-55 at pharmacologically relevant doses are primarily TopI poisons with DNA cleavage patterns exhibiting similarities and differences from those of CPT, but they are much more effective in *Leishmania* than in human enzyme.

342 **Induction of TopI-DNA complexes by indotecan and AM13-55 in *L. infantum***
343 **cultures**

344 To determine whether indotecan and AM13-55 induce TopI cleavage complexes in
345 drug-treated cells, we used the SDS-KCl precipitation assay. IFP1.4-*L. infantum*
346 (BCN150) promastigotes were exposed for 30 min to different concentrations of CPT
347 (used as positive control), indotecan and AM13-55 in a concentration range from 0.1 to
348 90 μ M. After this time, stabilized protein-DNA complexes were quantified by the
349 SDS/KCl precipitation method. Fig. 3A shows that CPT produced an increasing amount
350 of SDS/KCl-precipitable complexes that were dependent on the concentration of drug
351 that was added up to a value close to 80% of total labeled DNA. None of the
352 indenoisoquinolines studied had such potent effect within the same concentration range.
353 Indotecan and AM13-55 clearly showed an increase in the amount of SDS/KCl-
354 precipitable complexes, but unlike CPT, indotecan produced up to 20% labeled DNA at
355 10 μ M. AM13-55 reached up to 40% at the same concentration. These results are higher
356 than those obtained by Bakshi and co-workers with three sets of indenoisoquinolines
357 against trypanomastigotes of *T. brucei* (4). In that case, the indenoisoquinolines tested
358 only induced up to 12% of cleavage complexes from the total labeled DNA.

359 Since indotecan and AM13-55 inhibit DNA relaxation and induce stabilization of DNA
360 cleavage complexes, we performed a competition study between CPT and both
361 indenoisoquinolines in order to evaluate the primary mediator of cell killing in
362 Leishmania. Fig. 3B shows that 5 μ M of both tested compounds were not able to
363 prevent CPT-mediated TopIB-DNA stabilization thus pointing to a poison nature of
364 both drugs rather than interaction to DNA.

365 In order to assess the arresting effect of indotecan and AM13-55 on DNA synthesis, $5 \times$
366 10^6 exponentially growing *L. infantum* promastigotes were pulsed with 0.5 μ Ci [$2\text{-}^{14}\text{C}$]
367 thymidine in the presence of 1 μ M CPT, 1 μ M AM13-55, 1 μ M indotecan or solvent
368 (control). After a time lapse of 2, 4, 6, 8, 10 and 24 h, labeled DNA was determined by
369 scintillation counting. All the TopI poisons arrested DNA synthesis at all time points in
370 more than 90% with no significant differences amongst them (Fig. 3C). This results
371 correlate well with those found by Cushman *et al.* in 2011. They showed that
372 indenoisoquinolones with large primary amines side chain like indotecan and AM13-55 do
373 not intercalate into free DNA and do not suppress cleavage complex formation at high
374 concentration, unlike those with small primary amine (4,19).

375

376 **Indotecan and AM13-55 used as therapeutic agents in a visceral leishmaniasis**
377 **murine model infected with *L. infantum***

378 To evaluate the therapeutic potential of indotecan and AM13-55 *in vivo*, we performed
379 an experimental infection in a susceptible BALB/c strain mouse with *L. infantum*
380 promastigotes (wild type strain). Seven million infective-stage metacyclic promastigotes
381 were administered intravenously by the caudal vein of 15 healthy mice that were four to
382 six weeks old. By day 15 post-parasite inoculation, mice were divided into three groups
383 of five animals each. Mice were treated with 0.5 ml solutions of indotecan or AM13-55
384 in DMSO/saline, equivalent to 2.5 mg/kg body weight per injection, intraperitoneally
385 every two days for 15 days (total eight doses). As control, indenoisoquinoline vehicle
386 was administered.

387 Five days after the last administration, animals were killed by cervical dislocation and
388 spleens and livers were aseptically dissected to determine the parasite load. This was
389 determined by the limiting dilution assay. Each treatment group was compared to the
390 control, which had received only the vehicle in which the drug had been dissolved.

391 Fig. 4A shows the parasite burden in the group of mice treated with 2.5 mg/kg of
392 indotecan in spleen (upper panel) and liver (lower panel). After the administration
393 schedule was completed, a drastic reduction ($p<0.001$) of the number of transforming
394 amastigotes recovered from the target organs of drug-treated animals was observed. On
395 the other hand, mice treated with the same dose of AM13-55 (fig 4B) showed that only
396 spleen was efficiently cleared of infecting parasites, unlike liver, which retained the
397 pathogen load at similar levels to the untreated group (Fig 4B). In a subsequent trial, 5.0
398 mg/kg of body weight of AM13-55 was administered under the same conditions as
399 previous experiments (data not shown). Parasite load was reduced above 90% in the
400 spleen, but not in the liver, which remained unchanged. The resistance of liver
401 macrophages to kill the parasites may be due to metabolic transformations of the parent
402 compound in the hepatic parenchyma on TopIB-inactive byproducts.

403 Despite the fact that no definitive cure was achieved with indotecan, our results are very
404 similar to those found with 20 mg/kg body weight of paromomycin alone or in
405 combination with 200 mg/kg body weight glucantime on experimental infections of
406 BALB/c visceral leishmaniasis (14). Furthermore, both indotecan and AM13-55 were

407 more effective than CPT (2.5 mg/kg body weight), both free and liposome-
408 encapsulated, in a murine model of *L. donovani* leishmaniasis. The authors of that study
409 found that parasite load in livers and spleens were just reduced a 55% average when
410 infected mice were treated with this TopI poison (24).

411 In conclusion, we have proven that the two indenoisoquinolines analyzed have strong
412 leishmanicidal activity both *in vitro* (*ex vivo* splenic explant cultures) and *in vivo*
413 (visceral leishmaniosis murine model). These compounds act by stabilizing the DNA-
414 LiTopIB cleavage complexes and inhibiting the intrinsic relaxation activity of human
415 and leishmanial TopIB enzymes. Indotecan has a very high selectivity index with
416 respect to host cells and depleted the parasitic burden of spleen and liver. These facts
417 make this compound an exceptional candidate for its development as new
418 leishmanicidal drug with a better therapeutic profile than others currently in use.

419

420 **ACKNOWLEDGEMENTS**

421 This research was supported by Ministerio de Ciencia y Tecnología (grants
422 AGL2009-11935/GAN and AGL2010-16078/GAN), Instituto de Salud Carlos III (grant
423 PI09/0448 and the Network of Tropical Diseases RICET RD06/0021/1004) and Junta
424 de Castilla y León (grant Gr238). RAV, CFP and ECA are pre-doctoral fellows granted
425 by RICET (ISCIII), Junta de Castilla y León (ESF; European Social Founding) and
426 University of León, respectively. This work was also made possible by the National
427 Institutes of Health (NIH) through support with Research Grant U01 CA89566.

428

429

430

431

432

433

434

435

436 **CAPTIONS TO FIGURES**

437 **Figure 1.** Chemical structures of the indenoisoquinolines indotecan and AM13-55 used
438 in the present study.

439 **Figure 2. Indotecan and AM13-55 induce TopI-DNA covalent complexes and**
440 **inhibit supercoiled DNA relaxation.** A) Supercoiled plasmid DNA TopIB-mediated
441 cleavage assay showing the displacement towards cleavage complexes mediated by both
442 indenoisoquinolines and CPT on LiTopIB. Samples were run on a 1% agarose gel
443 containing ethidium bromide to a final concentration of 40 pg/ μ l in order to separate
444 supercoiled and relaxed DNA. The results are representative of three independent trials.
445 B) Inhibition of supercoiled DNA relaxation by indotecan (top gel) and AM13-55
446 (bottom gel) mediated by human (left lanes) and leishmanial TopIB (right lanes).
447 Reaction mixtures were incubated at 37 °C in a final concentration of 150 μ M KCl and
448 then stopped with SDS up to a final concentration of 1% of reaction volume. Products
449 were resolved in a 1% agarose gel and visualized by ethidium bromide staining. The
450 results are representative of three independent trials.

451 **Figure 3. Induction of TopI-DNA complexes by indotecan and AM13-55 in *L.***
452 ***infantum* cultures and DNA synthesis inhibition.** SDS/KCl-precipitable enzyme-
453 DNA complexes at increasing concentrations of the drugs under study: A) CPT,
454 indotecan and AM13-55, in promastigotes of *L. infantum* after 30 min growth in the
455 presence of different concentrations of drugs. Notice that in CPT the x-scale is up to 10
456 μ M, whereas in panels indotecan and AM13-55 is up to 90 μ M. Results are expressed as
457 mean \pm SE of at least three different experiments in duplicate; B) indotecan and AM13-
458 55 increase CPT-induced cleavable complex formation. Leishmania promastigotes were
459 treated with 5 μ M indotecan, 5 μ M AM13-55 or 10 μ M CPT for 30 min, or with 5 μ M
460 indotecan or AM13-55 for 5 min prior to addition of 10 μ M CPT, followed by an
461 additional 25-min incubation; (*) $P < 0.01$; (**) $P < 0.05$ using paired t-Student test; C) [2-
462 14 C] thymidine incorporation into DNA of growing promastigotes in the presence of 1
463 μ M CPT (empty dots), 1 μ M AM13-55 (empty triangles), 1 μ M indotecan (solid
464 triangles) or solvent (solid dots). The amount of labeled DNA is expressed as total cpm
465 per assay at different time points (2, 4, 6, 8, 10 and 24 h). Results are expressed as mean
466 \pm SE of at least three different experiments in duplicate.

467 **Figure 4.** Indenoisoquinolines cleared tissue parasitic burden in experimentally infected
468 BALB/c mice with *L. infantum*. Three groups of five animals each were challenged with
469 10^7 metacyclic promastigotes administered by the caudal vein. Fifteen days after
470 infection animals were injected intraperitoneally every two days for a total of fifteen
471 days (eight doses total), with a dose of 2.5-mg/kg body weight per injection of A)
472 indotecan, B) AM13-55 or the corresponding vehicle (a solution of DMSO in sterilized
473 saline solution). Mice were killed five days after the last treatment and the spleens and
474 livers were aseptically removed, weighted and homogenized in medium supplemented
475 with 20 % (v/v) FBS. After 10 days, transforming promastigotes were counted and the
476 limit dilution was considered for determining the parasitic burden of each organ. The
477 results are representative of two independent trials; statistical differences were observed
478 between groups (*) $P < 0.001$ using paired t-Student test.

479

480

481

482

483

484

485

486

487

488

489

490

491

492

493

494

495

496

497

498

REFERENCES

1. Aït-Oudhia K, Gazanion E, Vergnes B, Oury B, Sereno D. 2011. Leishmania antimony resistance: what we know what we can learn from the field. *Parasitol. Res.* 109: 1225-1232.
2. Antony S, Kohlhagen G, Agama K, Jayaraman M, Cao S, Durrani FA, Rustum YM, Cushman M, Pommier Y. 2005. Cellular topoisomerase I inhibition and antiproliferative activity by MJ-III-65 (NSC 706744), an indenoisoquinoline topoisomerase I poison. *Mol. Pharmacol.* 67:523-530.
3. Bakshi RP, Shapiro TA. 2003. DNA topoisomerases as targets for antiprotozoal therapy. *Mini Rev. Med. Chem.* 3: 597-608.
4. Bakshi RP, Sang D, Morrell A, Cushman M, Shapiro TA. 2009. Activity of indenoisoquinolines against African trypanosomes. *Antimicrob. Agents Chemother.* 53:123-128.
5. Balaña-Fouce R, Garcia-Estrada C, Perez-Pertejo Y, Reguera RM. 2008. Gene disruption of the DNA topoisomerase IB small subunit induces a non-viable phenotype in the hemoflagellate *Leishmania major*. *BMC Microbiol.* 8:113.
6. Balaña-Fouce R, Redondo CM, Pérez-Pertejo Y, Díaz-González R, Reguera RM. 2006. Targeting atypical trypanosomatid DNA topoisomerase I. *Drug Discov. Today* 11:733-740.
7. Balaña-Fouce R, Reguera RM, Cubría JC, Ordoñez D. 1998. The pharmacology of leishmaniasis. *Gen. Pharmacol.* 30: 435-443.
8. Bhattacharya SK, Sinha PK, Sundar S, Thakur CP, Jha TK, Pandey K, Das VR, Kumar N, Lal C, Verma N, Singh VP, Ranjan A, Verma RB, Anders G, Sindermann H, Ganguly NK. 2007. Phase 4 trial of miltefosine for the treatment of Indian visceral leishmaniasis. *J. Infect. Dis.* 196: 591-598.
9. Berman J. 2005. Miltefosine to treat leishmaniasis. *Expert Opin. Pharmacother.* 6: 1381-1388.
10. Bodley AL, Shapiro TA. 1995. Molecular and cytotoxic effects of camptothecin, a topoisomerase I inhibitor, on trypanosomes and *Leishmania*. *Proc. Natl. Acad. Sci. U.S.A.* 92:3726-3730.

- 530 11. Bodley AL, Chakraborty AK, Xie S, Burri C, Shapiro TA. 2003. An unusual type
531 IB topoisomerase from African trypanosomes. *Proc. Natl. Acad. Sci. U.S.A.*
532 100:7539-7544.
- 533 12. Díaz-González R, Pérez-Pertejo Y, Ordóñez D, Balaña-Fouce R, Reguera RM.
534 2007. Deletion study of DNA topoisomerase IB from *Leishmania donovani*:
535 searching for a minimal functional heterodimer. *PLoS One* 2:e1177.
- 536 13. Díaz-González R, Pérez-Pertejo Y, Pommier Y, Balaña-Fouce R, Reguera RM.
537 2008. Mutational study of the "catalytic tetrad" of DNA topoisomerase IB from
538 the hemoflagellate *Leishmania donovani*: Role of Asp-353 and Asn-221 in
539 camptothecin resistance. *Biochem. Pharmacol.* 76:608-619.
- 540 14. Gangneux JP, Sulahian A, Garin YJ, Derouin F. 1997. Efficacy of aminosidine
541 administered alone or in combination with meglumine antimoniate for the
542 treatment of experimental visceral leishmaniasis caused by *Leishmania infantum*.
543 *J. Antimicrob. Chemother.* 40:287-289.
- 544 15. Gietz RD, Schiestl RH. 1991. Applications of high efficiency lithium acetate
545 transformation of intact yeast cells using single-stranded nucleic acids as carrier.
546 *Yeast* 7:253-263.
- 547 16. Hotez PJ, Molyneux DH, Fenwick A, Ottesen E, Ehrlich Sachs S, Sachs JD. 2006.
548 Incorporating a rapid-impact package for neglected tropical diseases with
549 programs for HIV/AIDS, tuberculosis, and malaria. *PLoS Med.* 3:e102.
- 550 17. Hsiang YH, Hertzberg R, Hecht S, Liu LF. 1985. Camptothecin induces protein-
551 linked DNA break via mammalian DNA topoisomerase I. *J. Biol. Chem.*
552 260:14873–14878.
- 553 18. Kimutai A, Ngure K, Kiprotich Tonui W, Muita Gicheru M, Bonareri Nyamwamu
554 L. 2009. Leishmaniasis in Northern and Western Africa: a review. *Afr. J. Infect.*
555 *Dis.* 3:14 – 25.
- 556 19. Kiselev E, DeGuire S, Morrell A, Agama K, Dexheimer TS, Pommier Y,
557 Cushman M. 2011. 7-azaindonoisoquinolines as topoisomerase I inhibitors and
558 potential anticancer agents. *J Med Chem.* 54:6106-6116.

- 559 20. Miró G, Cardoso L, Pennisi MG, Oliva G, Baneth G. 2008. Canine leishmaniosis
560 – new concepts and insights on an expanding zoonosis: part two. Trends Parasitol.
561 24: 371-377.
- 562 21. Ossorio V, Travi BL, Rensio AR, Peniche AG, Melby PC. 2011. Identification of
563 small molecule lead compounds for visceral leishmaniasis using novel *ex vivo*
564 splenic explant model system. Plos Negl. Trop. Dis. 5: e962.
- 565 22. Pommier Y. 2006. Topoisomerase I inhibitors: camptothecins and beyond. Nat.
566 Rev. Cancer. 6:789-802.
- 567 23. Prada CF, Balaña-Fouce R, Álvarez-Velilla R, Calvo-Álvarez E, Requena JM,
568 Desideri A, Pérez-Pertejo Y, Reguera RM. 2012. Gimatecan and other
569 camptothecin derivatives poison Leishmania DNA-topoisomerase IB leading to a
570 strong leishmanicidal effect. J. Antimicrob. Chemother. (in press).
- 571 24. Proulx ME, Desormeaux A, Marquis JF, Olivier M, Bergeron MG. 2001.
572 Treatment of visceral leishmaniasis with sterically stabilized liposomes containing
573 camptothecin, Antimicrob. Agents Chemother. 45:2623–2627.
- 574 25. Prudhomme M. 2003. Rebeccamycin analogues as anti-cancer agents. Eur. J.
575 Med. Chem. 38:123-140.
- 576 26. Ready PD. 2010. Leishmaniasis emergence in Europe. Euro Surveill. 15:19505.
- 577 27. Robinson KA, Beverley SM. 2003. Improvements in transfection efficiency and
578 tests of RNA interference (RNAi) approaches in the protozoan parasite
579 *Leishmania*. Mol. Biochem. Parasitol. 128: 217–228.
- 580 28. Sacks D, Anderson C. 2004. Re-examination of the immunosuppressive
581 mechanisms mediating non-cure of Leishmania infection in mice. Immunol. Rev.
582 301: 225-238.
- 583 29. Sacks DL, Perkins PV. 1984. Identification of an infective stage of *Leishmania*
584 promastigotes. Science 223:1417-1419.
- 585 30. Schwartz E, Hatz C, Blum J. 2006. New world cutaneous leishmaniasis in
586 travelers. Lancet Infect. Dis. 6: 342-349.

- 587 31. Shu X, Royant A, Lin MZ, Aguilera TA, Lev-Ram V, Steinbach PA, Tsien RY.
588 2009. Mammalian expression of infrared fluorescent proteins engineered from a
589 bacterial phytochrome. *Science* 324: 804-807.
- 590 32. Teicher AB. 2008. Next generation topoisomerase I inhibitors: Rationale and
591 biomarker strategies. *Biochem. Pharmacol.* 75:1262-1271.
- 592 33. Thrash C, Voelkel K, DiNardo S, Sternglanz R. 1984. Identification of
593 *Saccharomyces cerevisiae* mutants deficient in DNA topoisomerase I activity. *J.*
594 *Biol. Chem.* 259: 1375-1377.
- 595 34. Titus RG, Marchand M, Boon T, Louis JA. 1985. A limiting dilution assay for
596 quantifying *Leishmania major* in tissues of infected mice. *Parasite Immunol.* 7:
597 545-555.
- 598 35. Villa H, Otero-Marcos AR, Reguera RM, Balaña-Fouce R, Garcia-Estrada C,
599 Pérez-Pertejo Y, Tekwani BL, Myler PJ, Stuart KD, Bjornsti MA, Ordóñez D.
600 2003. A novel active DNA topoisomerase I in *Leishmania donovani*. *J. Biol.*
601 *Chem.* 278: 3521-3526.
- 602 36. Zuma AA, Cavalcanti DP, Maia MC, de Souza W, Motta MC. 2011. Effect of
603 topoisomerase inhibitors and DNA-binding drugs on the cell proliferation and
604 ultrastructure of *Trypanosoma cruzi*. *Int. J. Antimicrob Agents.* 37:449-456.
- 605
606
607
608
609
610
611
612
613
614

615 **Table 1.** IC₅₀ calculation after a 48-h period of exposure to the selected compounds of
 616 *L. infantum* promastigotes, *ex vivo* infected splenocytes and uninfected splenocyte under
 617 methods described in the corresponding section. IC₅₀ at 48 h of the compounds selected
 618 in freshly uninfected splenocyte culture and values of SI_{48h} between this cell line and
 619 infected splenocytes with *L. infantum* amastigotes *ex vivo* were calculated from the
 620 dose-response curves performed in triplicate in separate experiments after performing a
 621 nonlinear fitting with the SigmaPlot® program.

622

drug	IC ₅₀ <i>L. infantum</i> promastigotes	IC ₅₀ amastigotes	IC ₅₀ uninfected splenocyte culture	SI _{48h} *
Paromomycin	42.41 μM ± 1.65	9.20 μM ± 0.01	15.70 mM ± 3.54	1706.5
camptothecin (22)	1.12 μM ± 0.13	0.03 μM ± 0.01	0.62 μM ± 0.13	20.7
indotecan	0.10 μM ± 0.08	0.10 μM ± 0.05	57.16 μM ± 6.01	571.6
AM13-55	1.02 μM ± 0.09	0.10 μM ± 0.37	48.37 μM ± 3.68	483.7

623

624 * selectivity index; SI_{48h} = IC₅₀ amastigotes / IC₅₀ infected splenocytes

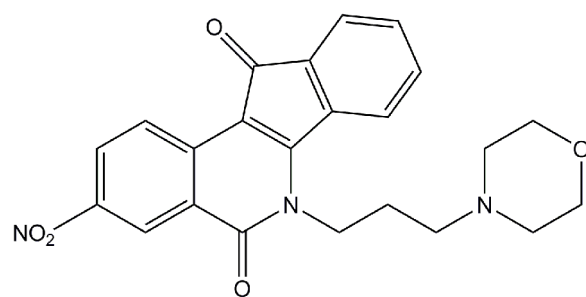
625

626

627

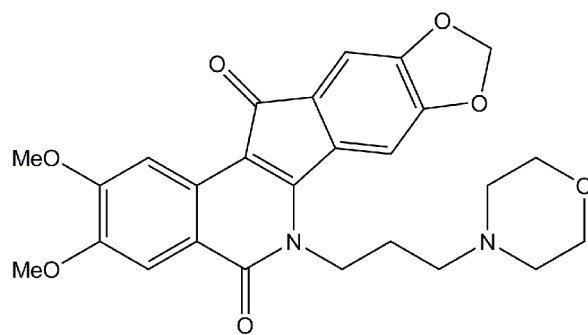
628

A



AM13-55

B



indotecan (LMP400)

