

Genome filtering identifies species-specific DNA biomarkers for *Mansonella perstans* and *Mansonella ozzardi* which enable differentiation of these closely related species and other co-endemic filarial parasites

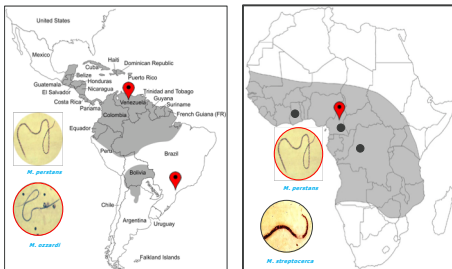
C. B. Poole¹, A. Sinha¹, L. Ettwiller¹, L. Apone¹, K. McKay¹, V. Panchapakesa¹, N. F. Lima², M. U. Ferreira², S. Wanji³ and C. K. S. Carlow¹

¹New England Biolabs, Ipswich, MA, USA. ²University of São Paulo, Brazil. ³University of Buea, Cameroon.

Introduction to Mansonelliasis

- Caused by 3 parasites: *Mansonella perstans*, *M. ozzardi* & *M. streptocerca*.
- Primary Insect Vector: *Culicoides* (biting midges).
 - *M. ozzardi* also spread by *Simulium* (black flies).
- No distinct, specific clinical consequences for *Mansonella* infections.
 - Immunosuppression caused by parasitic infection may lead to worsening of other medical conditions.
- Anti-helminthic treatment is complicated:
 - Not all species respond to ivermectin.
 - Benzimidazoles & DEC often employed.
- *Mansonella* patients are often co-infected with multiple filarial parasites including *Onchocerca volvulus*, *Wuchereria bancrofti*, and *Loa loa*.
- Need improved diagnostic tools.

Regions where Mansonelliasis is Endemic

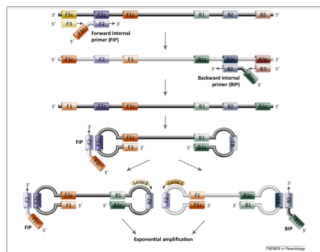


Grey shading represents regions endemic for Mansonelliasis

● Origin of parasite sample

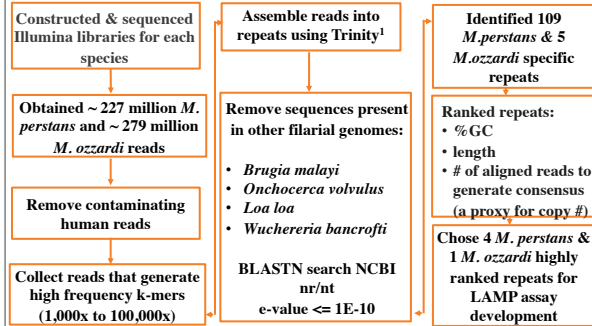
● Region co-endemic for *M. perstans* & *M. streptocerca*

Loop-Mediated Isothermal Amplification (LAMP) is Ideal for Field Use



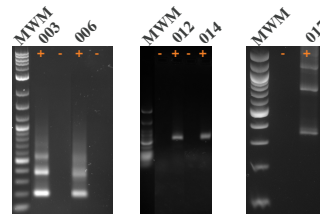
- Set of 4 primers (F3, FIP, B3 & BIP) recognize 6 distinct sequences on target DNA. Optional Loop primers speed up amplification.
- Single step isothermal reaction requiring a simple water bath.
- Highly sensitive and specific.
- Requires strand displacing polymerase (*Bst*).
- Rapid (~30-60 min) compared to PCR.
- Multiple direct methods for easy visualization of results.

Bioinformatic Pipeline Identifies New Diagnostic Biomarkers for *M. perstans* and *M. ozzardi*



¹Grabherr et al. (2011) *Nature Biotechnology* 29:644-654

Validation of Bioinformatically Identified Repeats by PCR



- PCR of candidate biomarkers from *M. perstans* (003, 006, 012 and 014) or *M. ozzardi* (017) DNA. Ladder-like arrays (003, 006 and 017) suggest repeats are organized tandemly in the genome whereas single bands (012, 014) suggest a dispersed organization. + = DNA; - = Non-template control; MWM = molecular weight marker.

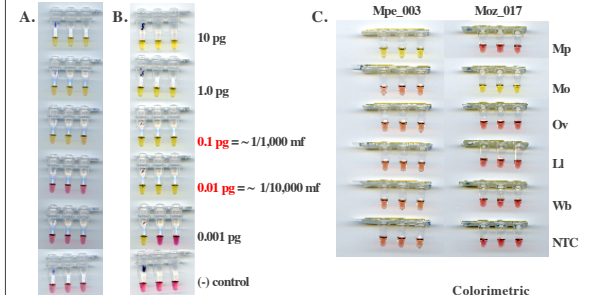
Candidate Biomarker Evaluation: Repeat Characteristics and LAMP Assay Results

Target	1/2 Reads aligned for consensus	%GC	Consensus sequence length (bps)	Repeat type	² Colorimetric LAMP		
					Optimal conditions	Sensitivity (pg)	Species specific
Mpe_003	80,315	33	366	tandem	63°C, 60 min	0.1	Yes
Mpe_006	111,796	31	579	tandem	63°C, 60 min	0.1	Yes
Mpe_012	21,974	29	338	dispersed	63°C, 60 min	1000	Yes
Mpe_014	13,223	29	318	dispersed	61°C, 60 min	10	Yes
Moz_017	10,000	42	303	tandem	63°C, 20 min	0.01	Yes

¹ # reads aligned to form the consensus sequence is used as a proxy for copy number.

² Multiple LAMP primer sets were designed for each target. The results generated by the best primer set are presented.

The *Mansonella* LAMP Assays are Sensitive and Specific



- The primer set targeting Mpe_003 detects as little as 0.1 pg *M. perstans* DNA (A) whereas the primer set targeting Moz_017 can detect as little as 0.01 pg of *M. ozzardi* DNA (B). The Mpe_003 and Moz_017 LAMP primer sets are specific for *M. perstans* (Mp) and *M. ozzardi* (Mo) DNA respectively. They do not cross react with DNA from the non-specific *Mansonella* species or with *O. volvulus* (Ov), *L. loa* (Ll) or *W. bancrofti* (Wb) DNA (C).

Validation on Patient and Insect Samples

Table 1. Detection of *M. perstans* in experimentally infected *C. milnei*. Comparison of the performance of ITS1 nested-PCR and colorimetric Mpe_003 LAMP.

<i>C. milnei</i> Infection Status	Sample Size	<i>Mansonella</i> Nested-PCR Positive	<i>M. perstans</i> LAMP Positive
FED ON VOLUNTEER: potentially infected	36	14	10
UNFED: presumed uninfected	36	1	0

Table 2. Detection of *M. perstans* in patient samples. Comparison of the performance of microscopy, ITS1 nested-PCR and colorimetric Mpe_003 LAMP.

Patient Infection Status ^a	Sample Size	<i>Mansonella</i> Nested-PCR Positive	<i>M. perstans</i> LAMP Positive
MF POSITIVE	9	1	9
MF NEGATIVE	1	1	1

^aAs determined by microscopic examination of patient samples.

Table 3. Detection of *M. ozzardi* in patient samples. Comparison of the performance of microscopy/ITS-2 qPCR, ITS1 nested-PCR and Moz_017 LAMP.

Patient Infection Status ^b	Sample Size	<i>Mansonella</i> Nested-PCR Positive	<i>M. ozzardi</i> LAMP Positive
MF / qPCR POSITIVE	51	Not Evaluated	51
MF / qPCR NEGATIVE	33	8	8

^bAs determined by microscopy and ITS-2 qPCR (Lima et al. (2018) *PLOS NTD* 12:e006327)

Summary

- Using a bioinformatic filtering approach, new diagnostic biomarkers for *M. perstans* and *M. ozzardi* were identified.
- Developed sensitive and species-specific LAMP assays targeting these new biomarkers.
- Validated these new LAMP assays on both patient and insect samples.
- These new field deployable assays will assist with the effort to better understand the global burden of Mansonelliasis.