# Malaria gametocytes production in the Wave Bioreactor: optimizing yields and quantification protocols



**ASTMH 2023 Meeting Poster Session A** 

Amel O.A. Ahmed, Standwell C. Nkhoma, Ioana Brasov, Sujatha Rashid, Rebecca Bradford and Robert E. Molestina American Type Culture Collection, 10801 University Boulevard, Manassas, VA 20110, USA Email: <u>aahmed@atcc.org</u>

Malaria is one of the most debilitating mosquito-borne diseases and poses a major health and socio-economic burden in many endemics of transmission are mainly determined by the density and sex ratio of the gametocytes. Molecular methods are critically needed for quantification of gametocytes, particularly when assessing transmission-blocking interventions. The lack of widely available and well-characterized gametocyte quantification reference standards hampers basic research on this crucial parasite stage and optimizing protocols for gametocyte quantification and sex discrimination. A known gametocyte-producing line MRA-1000 (NF-54), available from ATCC through BEI Resources, was used in optimization was performed using reverse transcriptase quantitative PCR assays targeting gametocyte-specific transcripts Pfs25 (female-specific), Pfs230 (male-specific), and Pfs16 (all gametocytes). Gametocytes). Gametocyte synthetic standards were generated by amplification of genomic DNA from a P. falciparum laboratory strain 3D7 (MRA-102) using sequence-specific primers containing a T7 promoter region that enables RNA transcription from in vitro DNA. RNA was transcribed into gametocytes can be accurately quantified and discriminated based on their sex using these synthetic doese in the sex using these synthetic contents and a discriminated based on their sex using these synthetic doese in the sex using these synthetic doese in the sex using the sex using these synthetic doese in the sex using terms are sex using the sex using terms are sex using te standards. These findings opened doors for research on this critical parasite stage. Availability of gametocyte-derived reagents and standards through BEI Resources.org), managed by ATCC, will aid research directed towards developing transmission-blocking interventions.

### BACKGROUND

**# LB-8021** 

- Transmission of *falciparum* malaria requires uptake of sexual stages (gametocytes) from infected individuals by mosquitoes. Gametocytes therefore pose as a logical target for malaria control and intervention studies.
- High-yield production of viable mature stage V gametocytes is a complex and time consuming process. Previously described protocols are labor and resource-intensive and usually generate limited quantities of gametocytes for research. This issue coupled with the absence of widely available and well-characterized gametocyte quantification standards hampers basic research on this critical parasite stage.
- The Wave Bioreactor presents an opportunity to produce vast quantities of gametocytes [1], in a controlled semi-automated system, which could be used for various applications including pre-clinical development of transmission blocking drugs and vaccines.
- The objectives of this study were to (i) optimize and establish a high-yield culture system for gametocyte production in the Wave 25 Bioreactor, and (ii) use synthetic in vitro cDNA standards developed in our lab to quantify and discriminate male from female gametocytes.

# **MATERIALS & METHODS**

#### In vitro culture of Plasmodium falciparum for gametocyte production

- Parasite strain. A known gametocyte-producing strain, NF-54 was obtained from the BEI Resources Repository, NIAID, NIH (<u>www.beiresources.org</u>): (BEI RESOURCES MRA-1000).
- **Parasite culture.** *P. falciprum* cultures were incubated at 37°C in leukocyte-depleted human type O+ erythrocytes (in CPDA-1) using RPMI 1640 supplemented with 10% heat-inactivated human type A+ serum, 0.18% Glucose, 0.18 mM Hypoxanthine, 1.77 mM L-Glutamine, 22 mM HEPES buffer 0.21% Sodium Bicarbonate and, 4 µg/mL Gentamicin under standard in vitro P. falciparum culture conditions
- Gametocyte induction and production. Gametocytes were produced in 2 L Cellbags on the Wave Bioreactor [1] (Figure 1). Induction of gametocytogenesis was performed in the cellbag upon inoculation at a high hematocrit (hct) and no media change for 48 hours.

#### **Detection and quantification of gametocyte-specific transcripts**

- Assays are based on detection of gametocyte-specific transcripts pfs16 (expressed in all gametocytes), pfs25 (female-specific) and pfs230 (male-specific) [2,3]
- Total RNA was extracted from gametocytemic cultures propagated in the Wave Bioreactor to make cDNA for RT-qPCR assays
- Synthetic in vitro DNA (ivDNA) standards were generated by amplifying genomic DNA from a standard *P. falciparum* laboratory strain 3D7 (BEI Resources www.beiresources.org) using sequence-specific primers containing a T7 promoter region that enables RNA transcription from in vitro DNA. RNA was transcribed into gametocyte-specific invitro cDNA (ivcDNA). The synthetic ivcDNA was used to generate standards for gametocyte quantification.
- RT-qPCR assays were performed using gametocyte-specific primers and probes (Life Technologies), and the CFX96<sup>™</sup> system (Bio-Rad Laboratories).

**AFS** 10801 University Boulevard, Manassas, VA 20110-2209

# ABSTRACT







#### **Phone:** 800.359.7370 **Email:** atccfederalsolutions@atcc.org

BEI Resources is funded by the National Institute of Allergy and Infectious Diseases (NIAID) and managed by ATCC. Materials are distributed only to established qualified research laboratories with facilities and safety programs appropriate for the level of material requested. © 2019 American Type Culture Collection. The ATCC trademark and trade name, and any other trademarks listed in this publication are trademarks owned by the American Type Culture Collection unless indicated otherwise





Figure 4. RT-qPCR Assay for detection and quantification of gametocytes. cDNA from harvested gametocytes was amplified and quantified using gametocyte-specific primers and probes. The assay is based on the detection of gametocyte-specific transcripts produced when parasites undergo sexual development (pfs16) – all gametocytes. Figures A and B show standard curves (generated from serially diluted ivcDNA synthetic standard for 3D7-pfs16. We extrapolated the quantity of gametocytes produced in a 2L Cellbag to be > 500 million gametocytes.



40

CellBag

Figure 5. Gametocyte quantification by microscopy.

Microscopy counts (100X magnification) of gametocytes in the WAVE bioreactor Cellbag (pelleted cells) and the supernatant (spent CB). Approximately 70% in gametocytes are females (green bars) and 30% males (blue bars). Gametocyte sex ratios tend to be female-biased.

We have optimized a protocol for high-yield production of malaria gametocytes in a semiautomated WAVE Bioreactor System.

• We developed and used synthetic gametocyte quantification standards to detect and quantify all gametocytes and discriminate the two gametocyte sexes. > 500 million gametocytes were harvested from the WAVE bioreactor.

These tools will help advance basic research on malaria gametocytes including understanding gametocyte sex ratio allocation and pre-clinical development of transmission-blocking drugs and vaccines.

Demanga C.G. et al. Parasit Vectors. 2017 May 2;10(1):216 2. Gadalla A.A.H. *et al*. Sci Rep. 2021 Sep 27;11(1):19118. 3. Schneider P. et al. Mol Biochem Parasitol. 2015 Jan-Feb;199(1-2):29-33

## ACKNOWLEDGEMENTS

This work was funded by the ATCC Internal Research and Development Program. The following strains were obtained from BEI Resources Repository, NIAID, NIH (www.beiresources.org): Plasmodium falciparum NF-54 (BEI

© ATCC 2019. The ATCC trademark, trade name, any and all ATCC catalog numbers listed in this publication are trademarks of the American Type Culture Collection unless indicated otherwise.