# DEI RESOURCES

# SUPPORTING INFECTIOUS DISEASE RESEARCH

ABSTRACT: Measurement of malaria parasite proliferation in cultured erythrocytes is critical for elucidating key determinants of phenotypes, including in complex growth interactions and variations in phenotype. We measured proliferation rates of three Plasmodium falciparum Cambodian isolates, including IPC\_3445 (MRA-1236), IPC\_5202 (MRA-1236), vitro cultures were maintained over four consecutive asexual parasite lineages, growth rates and relative changes in parasite mia were measured for laboratory parasite lines 3D7 (MRA-102) and DD2 (MRA-150) as controls. We observed significant differences in fold-change in parasite isolate MRA-1240 exhibits similar a proliferation rate to one of its constituent lineages, MRA1240-hap1 (GR: 1.03 ± 0.02 vs. 1.02 ± 0.05; FC: 67 ± 6 vs. 62 ± 2; p > 0.05), the other two component lineages (MRA1240-hap3) exhibit markedly different growth phenotype, masking the effect of minority haplotypes akin to recent observations from drug susceptibility testing. Our results also show diminished proliferation of isolate MRA-1236 (GR: 0.90 ± 0.02; FC: 101 ± 3) and MRA1236-hap2 (GR: 1.05 ± 0.04; FC: 88 ± 6) suggestive of competitive suppression. All parasite lines are available through BEI Resources and have well-defined in vitro growth phenotypes useful for both research and development of interventions against malaria.

### BACKGROUND

- Clinical malaria isolates often contain multiple parasite lineages exhibiting phenotypic variation
- Interaction amongst multiple lineages within a clinical isolate can drive evolution of important traits including growth rates, virulence and drug susceptibility
- \*We hypothesized that parasite lineages previously isolated from multiclonal isolates exhibit significantly variable proliferation rates under standard culture conditions
- To test this hypothesis, parasite growth rates and cell cycle times were measured for three Plasmodium falciparum Cambodian isolates, including IPC\_3445 (MRA-1236), IPC\_5202 (MRA-1240), IPC\_6403 (MRA-1285), and parasite lineages previously cloned from each of these isolates by limiting dilution<sup>1</sup>

# **MATERIALS & METHODS**

# *P. falciparum strains used and BEI Resources/MR4 catalog item numbers*

- P. falciparum IPC\_3445 (MRA-1236) and constituent haplotypes MRA1236-hap1 and MRA1236-hap2 ✤ P. falciparum IPC\_5202 (MRA-1240) and component haplotypes MRA1240-hap1, MRA1240-hap2, MRA1240hap3
- ✤ P. falciparum IPC 6403 (MRA-1285) and its three component haplotypes MRA1285-hap1, MRA1285-hap2 and MRA1285-hap3
- ✤ P. falciparum Dd2 (MRA-150)
- ✤ *P. falciparum* 3D7 (MRA-102)

# Standard conditions of asexual *P. falciparum* culture

All parasites were grown in leukocyte-depleted human type O+ erythrocytes at 37°C using a gas mixture containing 90% N<sub>2</sub>, 5% O<sub>2</sub> and 5% CO<sub>2</sub>

# **Composition of media used**

RPMI 1640 media (Gibco; Cat # 21870-084) supplemented with 4µg/mL Gentamicin solution (Gibco; Cat # 15750-060), 0.21% Sodium Bicarbonate (Gibco; Cat # 21870-084), 22mM HEPES buffer (Gibco; Cat # 15630-080), 0.18mM Hypoxanthine (Sigma; Cat # H9636), 0.18% Glucose (Sigma; Cat # G7021), 1.77mM L-Glutamine (Gibco; Cat # 25030-149) and 10% pooled human serum<sup>2</sup>

# Establishing highly synchronous cultures for growth rate studies

- ✤ Culture-adapted parasites (>3% parasitemia, >70% ring-stage) were synchronized by sorbitol treatment & incubated at 37°C for 42 hrs.
- \* Following incubation, cultured samples were passed through percoll gradients to enrich for schizonts
- Schizonts were grown for three hours & the resulting post-invasion rings (0-3 hr. old) were synchronized by sorbitol treatment, followed by multiple washes

### Setting up growth rate experiments

- ◆ 0-3 hr. old rings were seeded at 0.20% parasitemia and 3% hematocrit in 25cm<sup>2</sup> culture flasks & incubated at A 37°C for 48 hrs.
- Thin smears were made from each culture every 48 hrs. with fresh media changes made to each culture but without any blood added
- Smears were stained with 10% geimsa and read by examination of at least 500 cells with a light microscope

# Estimation of growth phenotypes

- ✤ For each parasite line we estimated both the fold-change in parasitemia & growth rate \* Fold-change describes change in parasitemia over time and is calculated as final parasitemia at the end of growth
- monitoring (in our case, after the 4<sup>th</sup> growth cycle) divided by the initial parasitemia used to seed the cultures Srowth rate per asexual cycle is estimated as the slope of a plot of log-transformed parasite counts against number of cycles

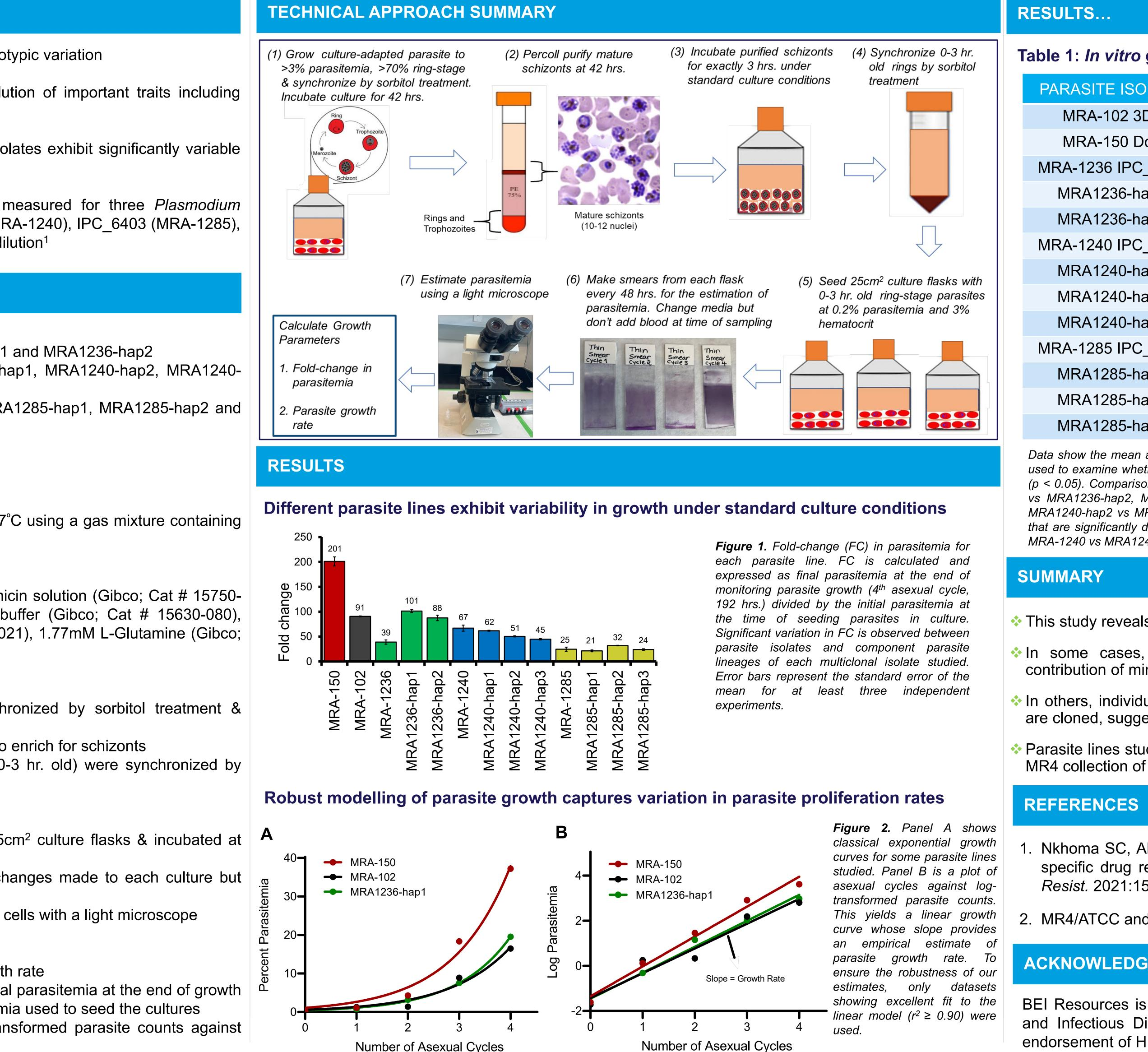




BEI Resources is funded by the National Institute of Allergy and Infectious Diseases (NIAID) and managed by ATCC. Materials are distributed only to established qualified research laboratories with facilities and safety programs appropriate for the level of material requested. © 2019 American Type Culture 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.

# In vitro growth phenotypes of single parasite lineages cloned from multiclonal malaria isolates

# Standwell C. Nkhoma, Amel O. Ahmed and Timothy T. Stedman BEI Resources, ATCC, 10801 University Boulevard, Manassas, VA 20110-2209, USA



# Number of Asexual Cycles

**Phone:** 800.359.7370

# 10801 University Boulevard, Manassas, Virginia 20110-2209

# **Email:** beiacquisitions@atcc.org

# **Web:** www.beiresources.org

### Table 1: In vitro growth phenotypes of parasite lines studied GROWTH RATE (MEAN ± SEM) PARASITE ISOLATE FOLD CHANGE (MEAN ± SEM) MRA-102 3D7 $1.08 \pm 0.01$ 91 ± 1 MRA-150 Dd2 201 ± 9 $1.27 \pm 0.04$ MRA-1236 IPC 3445 39 ± 4 $0.90 \pm 0.02$ 101 ± 3 $1.11 \pm 0.02$ MRA1236-hap1 MRA1236-hap2 88 ± 6 $1.01 \pm 0.04$ MRA-1240 IPC 5202 67 ± 6 $1.03 \pm 0.02$ 62 ± 2 MRA1240-hap1 $1.02 \pm 0.05$ MRA1240-hap2 51 ± 1 $0.94 \pm 0.01$ $0.93 \pm 0.01$ MRA1240-hap3 45 ± 1 $25 \pm 4$ $0.77 \pm 0.06$ MRA-1285 IPC 6403 MRA1285-hap1 21 ± 2 $0.80 \pm 0.02$ MRA1285-hap2 32 ± 1 $0.84 \pm 0.02$ MRA1285-hap3 24 ± 3 $0.82 \pm 0.04$

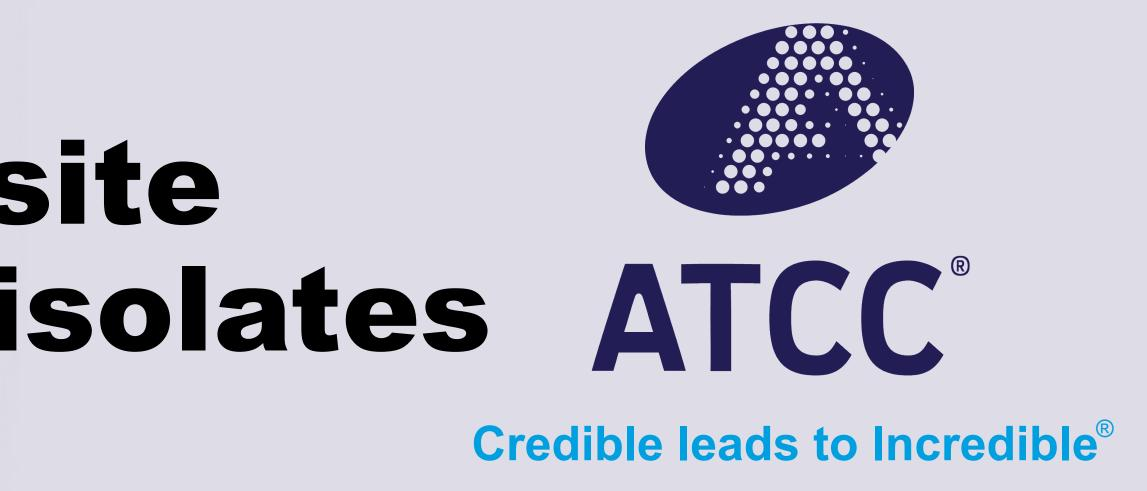
Data show the mean and standard error of the mean (SEM) for at least three independent experiments. An unpaired t-test was used to examine whether empirical values for the fold-change or growth rates of any two parasite lines are significantly different (p < 0.05). Comparisons that are significantly different for the mean fold-change are: MRA-1236 vs MRA1236-hap1, MRA-1236 vs MRA1236-hap2, MRA-1240 vs MRA1240-hap3, MRA1240-hap1 vs MRA1240-hap2, MRA1240-hap1 vs MRA1240-hap3, MRA1240-hap2 vs MRA1240-hap3, MRA1285-hap1 vs MRA1285-hap2 and MRA1285-hap2 vs MRA1285-hap3. Comparisons that are significantly different (p < 0.05) for the growth rate are: MRA-1236 vs MRA1236-hap1, MRA-1236 vs MRA1236-hap2, MRA-1240 vs MRA1240-hap2 and MRA-1240 vs MRA1240-hap3.

- contribution of minor frequency lineages

- Resist. 2021:152-161.

# ACKNOWLEDGEMENTS

BEI Resources is funded under contract HHSN272201600013C by the National Institute of Allergy and Infectious Diseases, NIH. The views expressed in this work do not necessarily reflect the endorsement of HHS.



This study reveals significant variation in proliferation between co-infecting parasite lineages

In some cases, the most predominant lineage dominates the growth profile, masking the

In others, individual parasite lineages exhibit greater proliferation than the isolate from which they are cloned, suggestive of competitive suppression

\* Parasite lines studied here are available to the malaria research and control community through the MR4 collection of NIAID's BEI Resources Program (https://www.beiresources.org)

Nkhoma SC, Ahmed AO, Zaman S, Porier D, Baker Z and Stedman TT. Dissection of haplotypespecific drug response phenotypes in multiclonal malaria isolates. Int. J. Parasitol. Drugs Drug

2. MR4/ATCC and EVIMalaR. Methods in Malaria Research. (2013) 6th ed., Manassas, VA.