
Histidine-Rich Protein PFHRP-II Polyclonal Antibody

(Catalog # A64450)

Description

Histidine-Rich Protein PFHRP-II Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

Specificity

Plasmodium falciparum

Isotype

IgG

Uniprot ID

P05227

Purification

>95%, Protein G purified

Immunogen

Recombinant Plasmodium falciparum Histidine-rich protein PFHRP-II protein (31-289AA)

Storage

Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Histidine-rich protein PFHRP-II

Application

ELISA
