
AMA-1 Polyclonal Antibody

(Catalog #A53015)

Background

Involved in parasite invasion of erythrocytes.

Description

AMA-1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

0.1mol/L NaCl, 15mmol/L NaN₃, PH7.2 .

Specificity

Plasmodium falciparum

Isotype

IgG

Uniprot ID

P22621

Purification

Caprylic Acid Ammonium Sulfate Precipitation purified

Immunogen

Recombinant Plasmodium falciparum Apical membrane antigen 1 protein (25-546AA)

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

Merozoite surface antigen, AMA-1

Application

ELISA
