
PRAMEF5 Polyclonal Antibody

(Catalog #A63214)

Description

PRAMEF5 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

Q5TYX0

Purification

>95%, Protein G purified

Immunogen

Recombinant Human PRAME family member 5 protein (1-476AA)

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

PRAME family member 5, PRAMEF5

Application

ELISA; Recommended dilution: ELISA: 1:500-1:1000
