
WFDC9 Polyclonal Antibody

(Catalog # A66402)

Description

WFDC9 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

Q8NEX5

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Protein WFDC9 protein (24-89AA)

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

Protein WFDC9, WFDC9, WAP9

Application

ELISA: Recommended dilution: 1:500 -1:2000
