

WFDC11 Polyclonal Antibody

(Catalog # A66222)

Description

WFDC11 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

Q8NEX6

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Protein WFDC11 protein (26-87AA)

Storage

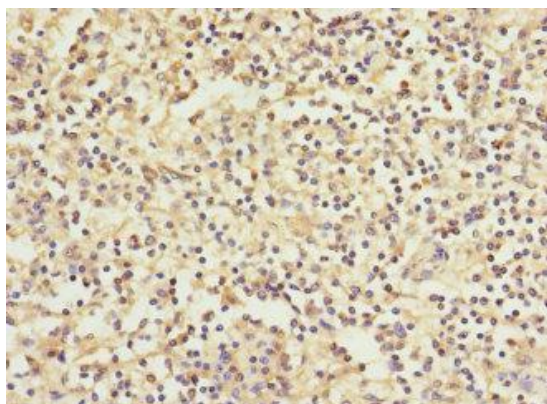
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Protein WFDC11, WFDC11, WAP11

Application

ELISA, IHC; Recommended dilution: IHC:1:20-1:200



Immunohistochemistry of paraffin-embedded human spleen tissue WFDC11 Polyclonal Antibody at dilution 1:100