
C22orf23 Polyclonal Antibody

(Catalog #A67462)

Description

C22orf23 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

Q9BZE7

Purification

>95%, Protein G purified

Immunogen

Recombinant Human UPF0193 protein EVG1 protein (1-217AA)

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

UPF0193 protein EVG1, C22orf23

Application

ELISA; Recommended dilution: ELISA: 1:1000
