
ALG14 Polyclonal Antibody, FITC Conjugated

(Catalog #A58052)

Background

May be involved in protein N-glycosylation. May play a role in the second step of the dolichol-linked oligosaccharide pathway. May anchor the catalytic subunit ALG13 to the ER.

Description

ALG14 Polyclonal Antibody, FITC Conjugated. FITC. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

Q96F25

Purification

>95%, Protein G purified

Immunogen

Recombinant Human UDP-N-acetylglucosamine transferase subunit ALG14 homolog protein (25-216AA)

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

UDP-N-acetylglucosamine transferase subunit ALG14 homolog, ALG14

Application

ELISA; Recommended dilution: ELISA 1:100-1:500
