

NACA2 Polyclonal Antibody

(Catalog # A67934)

Background

Prevents inappropriate targeting of non-secretory polypeptides to the endoplasmic reticulum (ER). Binds to nascent polypeptide chains as they emerge from the ribosome and blocks their interaction with the signal recognition particle (SRP), which normally targets nascent secretory peptides to the ER. Also reduces the inherent affinity of ribosomes for protein translocation sites in the ER membrane (M sites).

Description

NACA2 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

Q9H009

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Nascent polypeptide-associated complex subunit alpha-2 protein (1-215AA)

Storage

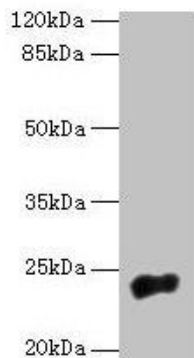
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Nascent polypeptide-associated complex subunit alpha-2, Alpha-NAC-like, Hom s 2.01, Nascent polypeptide-associated complex subunit alpha-like, NAC-alpha-like, NACA2, NACAL

Application

ELISA, WB; Recommended dilution: WB:1:1000-1:5000



Western blot

All lanes: NACA2 Polyclonal Antibody at 5ug/ml+293T whole cell lysate

Goat polyclonal to rabbit at 1/10000 dilution

Predicted band size: 23 kDa

Observed band size: 23 kDa