
NUDCD2 Polyclonal Antibody

(Catalog # A66662)

Background

May regulate the LIS1/dynein pathway by stabilizing LIS1 with Hsp90 chaperone.

Description

NUDCD2 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

Q8WVJ2

Purification

>95%, Protein G purified

Immunogen

Recombinant Human NudC domain-containing protein 2 protein (1-157AA)

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

NudC domain-containing protein 2, NUDCD2

Application

ELISA
