

DYNLRB1 Polyclonal Antibody

(Catalog # A51253)

Background

Acts as one of several non-catalytic accessory components of the cytoplasmic dynein 1 complex that are thought to be involved in linking dynein to cargos and to adapter proteins that regulate dynein function. Cytoplasmic dynein 1 acts as a motor for the intracellular retrograde motility of vesicles and organelles along microtubules.

Description

DYNLRB1 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

Q9NP97

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Dynein light chain roadblock-type 1 protein (3-96AA)

Storage

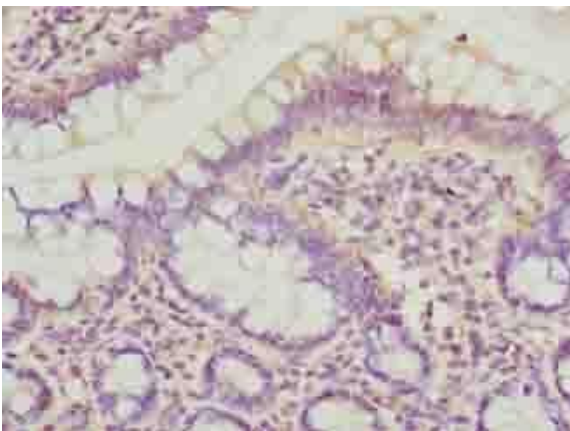
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

BLP, Dynein light chain 2A, cytoplasmic, Dynein-associated protein Km23, Roadblock domain-containing protein 1, BITH, DNCL2A, DNLC2A, ROBLD1, HSPC162, DYNLRB1

Application

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



Immunohistochemistry of paraffin-embedded human small intestine tissue using DYNLRB1 Polyclonal Antibody at dilution of 1:100