

TRAK1 Polyclonal Antibody

(Catalog A68042)

Background

Involved in the regulation of endosome-to-lysosome trafficking, including endocytic trafficking of EGF-EGFR complexes and GABA-A receptors.

Description

TRAK1 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

Q9UPV9

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Trafficking kinesin-binding protein 1 protein (101-300AA)

Storage

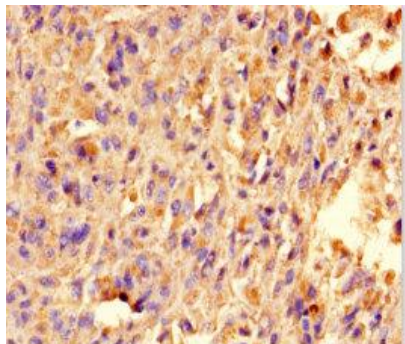
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Trafficking kinesin-binding protein 1, 106 kDa O-GlcNAc transferase-interacting protein, TRAK1, KIAA1042, OIP106

Application

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



Immunohistochemistry of paraffin-embedded human melanoma cancer using TRAK1 Polyclonal Antibody at dilution 1:100