
RHEBL1 Polyclonal Antibody

(Catalog #A60602)

Background

Binds GTP and exhibits intrinsic GTPase activity. May activate NF-kappa-B-mediated gene transcription. Promotes signal transduction through MTOR, activates RPS6KB1, and is a downstream target of the small GTPase-activating proteins TSC1 and TSC2.

Description

RHEBL1 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

Q8TAI7

Purification

>95%, Protein G purified

Immunogen

Recombinant Human GTPase RhebL1 protein (1-180AA)

Storage

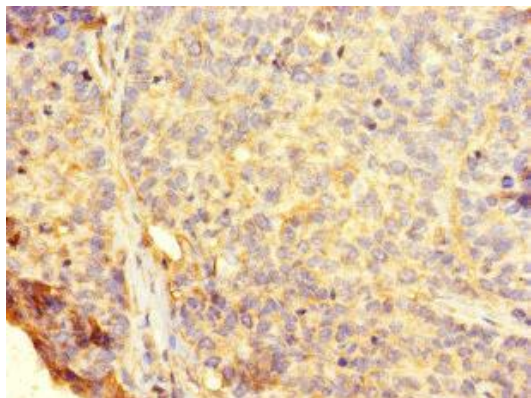
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

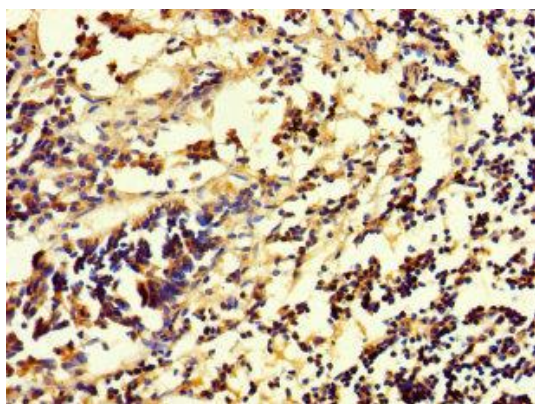
GTPase RhebL1, RHEBL1, Ras homolog enriched in brain like-1 c, RhebL1c, Ras homolog enriched in brain-like protein 1, Rheb-like protein 1, Rheb2

Application

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



Immunohistochemistry of paraffin-embedded human ovarian cancer using RHEBL1 Polyclonal Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human appendix tissue using RHEBL1 Polyclonal Antibody at dilution of 1:100