

TTLL6 Polyclonal Antibody

(Catalog # A66198)

Background

Polyglutamylase which preferentially modifies alpha-tubulin. Mediates tubulin polyglutamylation in cilia. Involved in the side-chain elongation step of the polyglutamylation reaction rather than in the initiation step.

Description

TTLL6 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

Q8N841

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Tubulin polyglutamylase TTLL6 protein (121-310AA)

Storage

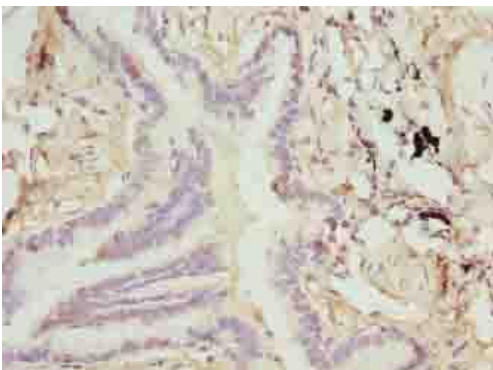
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Tubulin polyglutamylase TTLL6, Tubulin--tyrosine ligase-like protein 6, TTLL6

Application

ELISA, WB, IHC; Recommended dilution: WB:1:500-1:5000, IHC:1:500-1:1000



Immunohistochemistry of paraffin-embedded human lung tissue using TTLL6 Polyclonal Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human testis tissue using TTLL6 Polyclonal Antibody at dilution of 1:100