
TUBA1A Polyclonal Antibody, FITC Conjugated

(Catalog # A61612)

Background

Tubulin is the major constituent of microtubules. It binds two moles of GTP, one at an exchangeable site on the beta chain and one at a non-exchangeable site on the alpha chain.

Description

TUBA1A Polyclonal Antibody, FITC Conjugated. FITC. Raised in: Rabbit.

Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

Specificity

Human

Isotype

IgG

Uniprot ID

Q71U36

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Tubulin alpha-1A chain protein (30-229AA)

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 alpha-1A chain, TUBA1A, TUBA3, Alpha-tubulin 3, Tubulin B-alpha-1, Tubulin alpha-3 chain

Application

ELISA; Recommended dilution: ELISA 1:100-1:500
