
SPIN2A Polyclonal Antibody

(Catalog # A67146)

Background

May be involved in the regulation of cell cycle progression.

Description

SPIN2A 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

Q99865

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Spindlin-2A protein (1-258AA)

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

Spindlin-2A, Protein DXF34, Spindlin-like protein 2A, SPIN-2, SPIN-2A, SPIN2A, DXF34, SPIN2

Application

ELISA
