

SPIN3 Polyclonal Antibody

(Catalog # A65134)

Background

SPIN3 (Spindlin Family Member 3) is a Protein Coding gene that exhibits H3K4me3-binding activity. Diseases associated with SPIN3 include Atrial Heart Septal Defect. An important paralog of this gene is SPIN1.

Description

SPIN3 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

Q5JUX0

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Spindlin-3 protein (1-77AA)

Storage

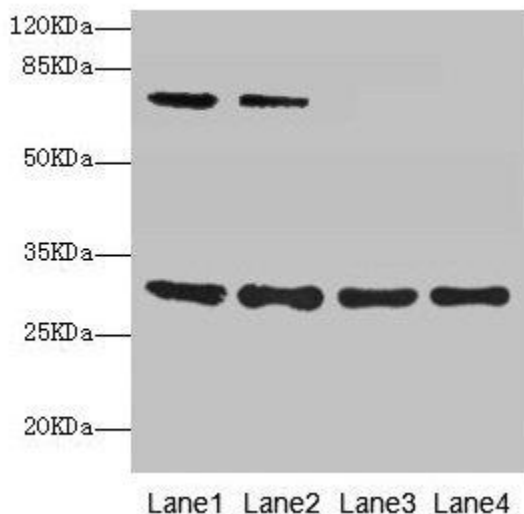
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Spindlin-3, Spindlin-like protein 3, SPIN-3, SPIN3

Application

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



Western blot All lanes: SPIN3 Polyclonal Antibody at 6µg/ml
Lane 1: A549 whole cell lysate
Lane 2: Hela whole cell lysate
Lane 3: A431 whole cell lysate
Lane 4: HepG2 whole cell lysate
Secondary
Goat polyclonal to rabbit IgG at 1/10000 dilution
Predicted band size: 30, 9 kDa
Observed band size: 30, 70 kDa