
SRP54 Polyclonal Antibody

(Catalog # A63498)

Background

Binds to the signal sequence of presecretory protein when they emerge from the ribosomes and transfers them to TRAM (translocating chain-associating membrane protein).

Description

SRP54 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

P61011

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Signal recognition particle 54 kDa protein (1-504AA)

Storage

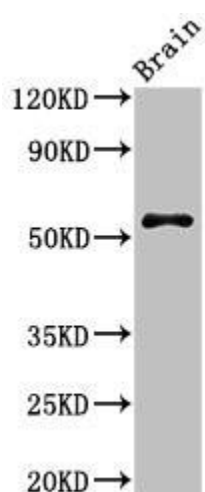
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Signal recognition particle 54 kDa protein, SRP54, SRP54

Application

ELISA, WB, IF; Recommended dilution: WB:1:500-1:5000, IF:1:50-1:200



Western Blot

Positive WB detected in: Mouse brain tissue

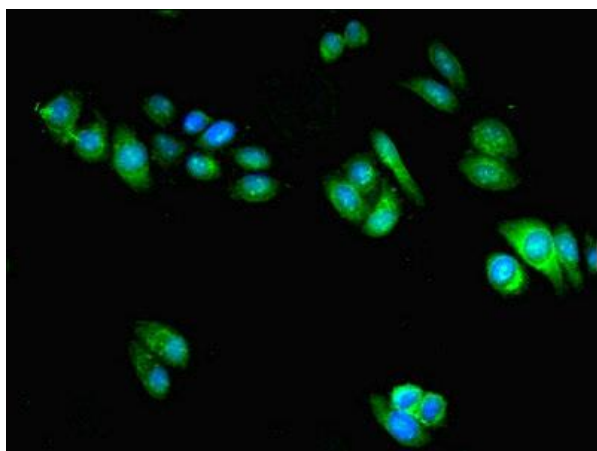
All lanes: SRP54 Polyclonal Antibody at 3ug/ml

Secondary

Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 56, 51 KDa

Observed band size: 56 KDa



Immunofluorescent analysis of HepG2 cells using SRP54

Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-
conjugated AffiniPure Goat Anti-Rabbit IgG (H+L).