

C10orf54 Polyclonal Antibody

(Catalog # A67834)

Background

May stimulate MMP14-mediated MMP2 activation.

Description

C10orf54 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

Q9H7M9

Purification

Protein G purified

Immunogen

Recombinant Human V-type immunoglobulin domain-containing suppressor of T-cell activation protein (37-194AA)

Storage

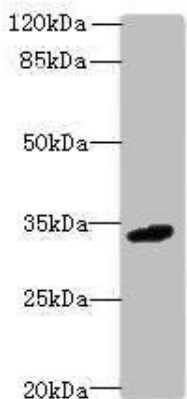
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Platelet receptor Gi24, Stress-induced secreted protein-1, Sisp-1, C10orf54, SISP1, PP2135, UNQ730/PRO1412, VSIR (V-set immunoregulatory receptor), VSIR

Application

ELISA, WB, IHC; Recommended dilution: WB:1:1000-1:5000, IHC:1:20-1:200



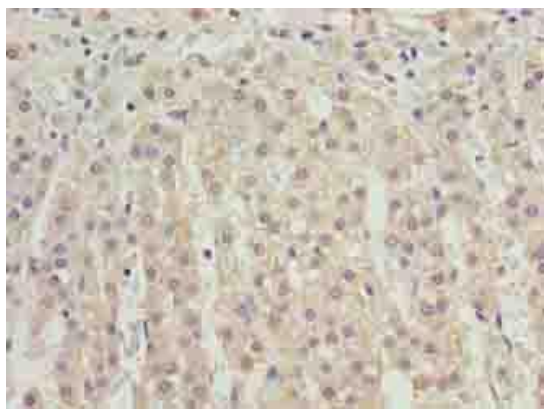
Western blot

All lanes: C10orf54 Polyclonal Antibody at 4ug/ml+Mouse liver tissue

Goat polyclonal to rabbit at 1/10000 dilution

Predicted band size: 34 kDa

Observed band size: 34 kDa



Immunohistochemistry of paraffin-embedded human adrenal gland tissue using C10orf54 Polyclonal Antibody at dilution of 1:100