

LRRC57 Polyclonal Antibody

(Catalog # A66762)

Background

extracellular exosome

Description

LRRC57 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

Q8N9N7

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Leucine-rich repeat-containing protein 57 protein (1-239AA)

Storage

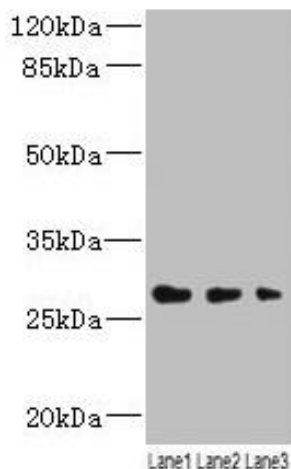
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Leucine-rich repeat-containing protein 57, LRRC57

Application

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



Western blot

All lanes: LRRC57 Polyclonal Antibody at 4 ug/ml

Lane 1: Mouse liver tissue

Lane 2: Mouse kidney tissue

Lane 3: Mouse lung tissue

Secondary

Goat polyclonal to Rabbit IgG at 1/10000 dilution

Predicted band size: 27 kDa

Observed band size: 27 kDa