

LRRC40 Polyclonal Antibody

(Catalog # A53788)

Description

LRRC40 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

Q9H9A6

Purification

Protein G purified

Immunogen

Recombinant Human Leucine-rich repeat-containing protein 40 protein (1-602AA)

Storage

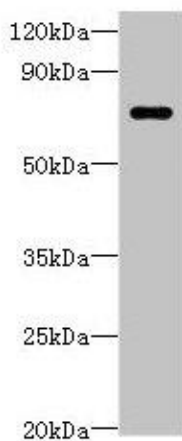
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

LRRC40

Application

ELISA, WB; Recommended dilution: WB:1:500-1:5000



Western blot

All lanes: LRRC40 Polyclonal Antibody at 2µg/ml + 293T whole cell lysate

Secondary

Goat polyclonal to rabbit IgG at 1/10000 dilution

Predicted band size: 69 kDa

Observed band size: 69 kDa