

LGALS12 Polyclonal Antibody

(Catalog # A59614)

Background

Binds lactose. May participate in the apoptosis of adipocytes.

Description

LGALS12 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Rat

Isotype

IgG

Uniprot ID

Q96DT0

Purification

Protein G purified

Immunogen

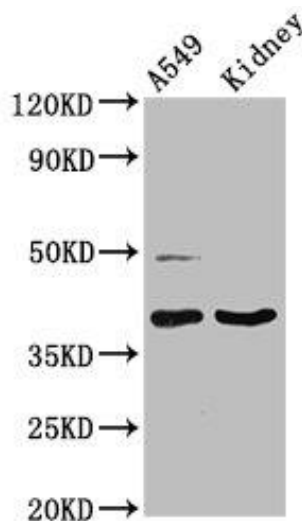
Recombinant Human Galectin-12 protein (1-336AA)

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

Galectin-12, Gal-12, LGALS12, GRIP1, Galectin-related inhibitor of proliferation



Western Blot

Positive WB detected in: A549 whole cell lysate, Rat kidney tissue

All lanes: LGALS12 Polyclonal Antibody at 3µg/ml
Secondary

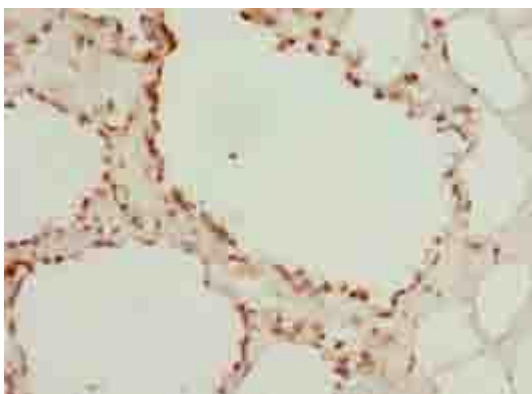
Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 38, 32, 37, 30, 36 kDa

Observed band size: 38 kDa



Immunohistochemistry of paraffin-embedded human skin tissue using LGALS12 Polyclonal Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human breast cancer using LGALS12 Polyclonal Antibody at dilution of 1:100