

FAM63A Polyclonal Antibody

(Catalog # A66742)

Background

extracellular exosome, nucleoplasm, nucleus

Description

FAM63A 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

Q8N5J2

Purification

Protein G purified

Immunogen

Recombinant Human Ubiquitin carboxyl-terminal hydrolase MINDY-1 protein (3-278AA)

Storage

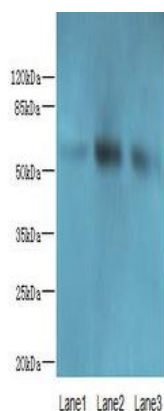
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Protein FAM63A, FAM63A, KIAA1390

Application

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



Western blot

All lanes: FAM63A Polyclonal Antibody at 8ug/ml

Lane 1: Mouse liver tissue

Lane 2: A549 whole cell lysate

Lane 3: HepG-2 whole cell lysate

Goat polyclonal to Rabbit IgG at 1/10000 dilution

Predicted band size: 52 kDa

Observed band size: 52 kDa