

KRT79 Polyclonal Antibody

(Catalog # A65214)

Background

extracellular exosome, enzyme binding

Description

KRT79 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

Q5XKE5

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Keratin, type II cytoskeletal 79 protein (1-300AA)

Storage

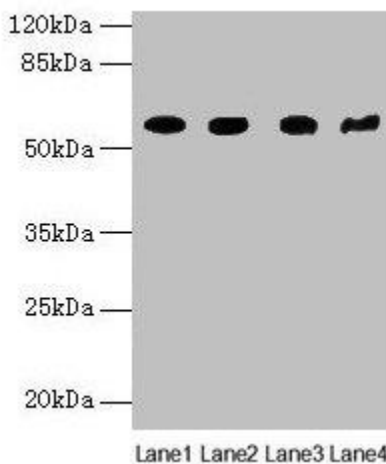
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Keratin, type II cytoskeletal 79, Cytokeratin-79, CK-79, Keratin-6-like, Keratin-6L, Keratin-79, K79, Type-II keratin Kb38, KRT79, K6L, KB38, KRT6L

Application

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



Western blot

All lanes: KRT79 Polyclonal Antibody at 1ug/ml

Lane 1: Mouse lung tissue

Lane 2: A549 whole cell lysate

Lane 3: 293T whole cell lysate

Lane 4: Mouse gonadal tissue

Goat polyclonal to Rabbit IgG at 1/10000 dilution

Predicted band size: 58 kDa

Observed band size: 58 kDa