

CTSL Polyclonal Antibody

(Catalog # A50161)

Background

Important for the overall degradation of proteins in lysosomes.

Description

CTSL Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Liquid. 0.03% Proclin 300. 50% Glycerol, 0.01M PBS, pH 7.4.

Specificity

Chlorocebus aethiops, Human

Isotype

IgG

Uniprot ID

Q9GKL8

Purification

Protein G purified

Immunogen

Recombinant Chlorocebus aethiops Cathepsin L1 protein (114-333AA)

Storage

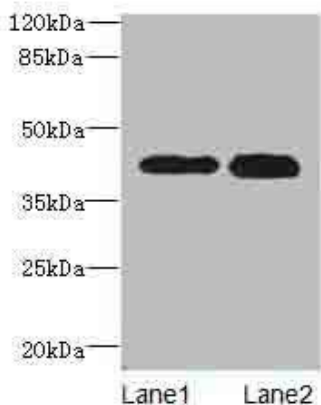
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Cathepsin L, Major excreted protein, MEP, CTSL1, CTSL

Application

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



Western blot

All lanes: CTSL Polyclonal Antibody at 12ug/ml

Lane 1: A549 whole cell lysate

Lane 2: HepG-2 whole cell lysate

Secondary

Goat polyclonal to Rabbit IgG at 1/10000 dilution

Predicted band size: 37 kDa

Observed band size: 37 kDa