

## ENTPD3 Polyclonal Antibody

(Catalog # A71977)

### Background

This gene encodes a plasma membrane-bound divalent cation-dependent E-type nucleotidase. The encoded protein is involved in the regulation of extracellular levels of ATP by hydrolysis of it and other nucleotides. Multiple transcript variants have been described.

### Description

ENTPD3 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

### Specificity

Mouse

### Isotype

IgG

### Uniprot ID

O75355

### Purification

Affinity Purification

### Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 45-366 of human ENTPD3 (NP\_001278890.1).

### Storage

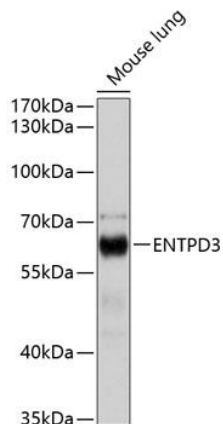
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

### Alternative Names

ENTPD3; CD39L3; HB6; NTPDase-3; ectonucleoside triphosphate diphosphohydrolase 3

### Application

WB; Recommended dilution: WB: 1:500-1:2000



#### Western Blot

Western blot analysis of extracts of mouse lung, using ENTPD3 Polyclonal Antibody at 1:1000 dilution.

Secondary Antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.

Lysates/proteins: 25ug per lane.

Blocking buffer: 3% nonfat dry milk in TBST.

Exposure time: 10s.