

## MCAM Polyclonal Antibody

(Catalog # A71130)

### Description

MCAM Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

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

### Specificity

Human, Rat

### Isotype

IgG

### Uniprot ID

P43121

### Purification

Affinity Purification

### Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 340-560 of human MCAM (NP\_006491.2).

### Storage

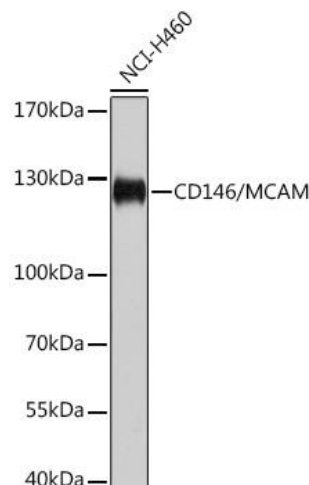
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

### Alternative Names

MCAM; CD146; MUC18; cell surface glycoprotein MUC18

### Application

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



Western blot analysis of extracts of NCI-H460 cells, using MCAM 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.