

## FERMT2 Polyclonal Antibody

(Catalog # A73197)

### Description

FERMT2 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

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

### Specificity

Human, Mouse, Rat

### Isotype

IgG

### Uniprot ID

Q96AC1

### Purification

Affinity Purification

### Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 380-550 of human FERMT2 (NP\_006823.1).

### Storage

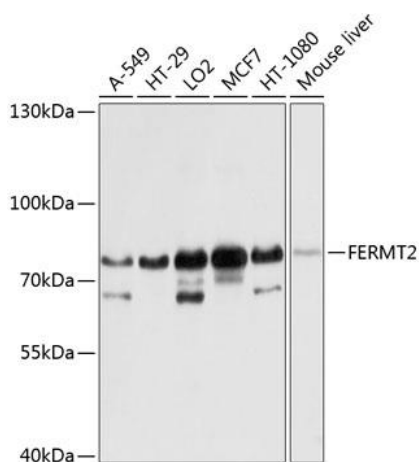
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

### Alternative Names

FERMT2; KIND2; MIG2; PLEKHC1; UNC112; UNC112B; mig-2; fermitin family member 2

### Application

WB, IHC; Recommended dilution: WB 1:1000 - 1:2000, IHC 1:50 - 1:200



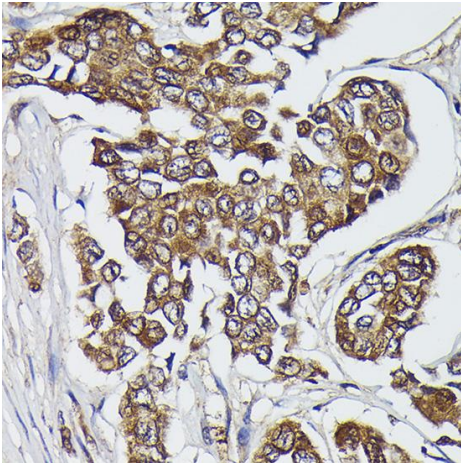
Western blot analysis of extracts of various cell lines, using FERMT2 Polyclonal Antibody at 1:1000 dilution.

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

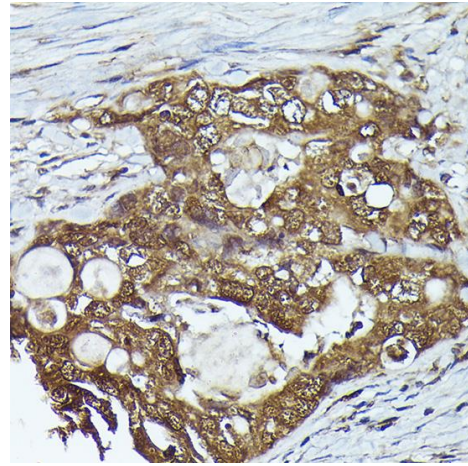
Lysates/proteins: 25ug per lane.

Blocking buffer: 3% nonfat dry milk in TBST.

Exposure time: 5s.



Immunohistochemistry analysis of paraffin-embedded human breast cancer using FERMT2 Rabbit pAb at dilution of 1:300 (40x lens). Performed high pressure antigen retrieval with 10 mM citrate buffer pH 6.0 before commencing with IHC staining protocol.



Immunohistochemistry analysis of paraffin-embedded human colon carcinoma using FERMT2 Rabbit pAb at dilution of 1:300 (40x lens). Performed high pressure antigen retrieval with 10 mM citrate buffer pH 6.0 before commencing with IHC staining protocol.