

## HM13 Polyclonal Antibody

(Catalog # A59386)

### Background

Catalyzes intramembrane proteolysis of some signal peptides after they have been cleaved from a preprotein, resulting in the release of the fragment from the ER membrane into the cytoplasm. Required to generate lymphocyte cell surface (HLA-E) epitopes derived from MHC class I signal peptides (PubMed:11714810). May be necessary for the removal of the signal peptide that remains attached to the hepatitis C virus core protein after the initial proteolytic processing of the polyprotein (PubMed:12145199). Involved in the intramembrane cleavage of the integral membrane protein PSEN1 (PubMed:12077416, PubMed:11714810, PubMed:14741365). Cleaves the integral membrane protein XBP1 isoform 1 in a DERL1/RNF139-dependent manner (PubMed:25239945). May play a role in graft rejection.

### Description

HM13 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

Q8TCT9

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human Minor histocompatibility antigen H13 protein (336-377AA)

### Storage

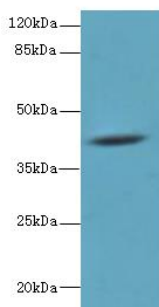
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Alternative Names

Minor histocompatibility antigen H13 (EC:3.4.23.-), HM13, H13, IMP1, PSL3, SPP, MSTP086, Intramembrane protease 1, IMP-1, IMPAS-1, hIMP1, Presenilin-like protein 3, Signal peptide peptidase

### Application

ELISA, WB, IF; Recommended dilution: WB:1:1000-1:5000, IF:1:50-1:200



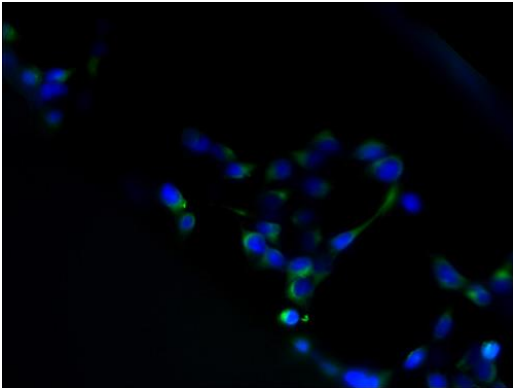
### Western blot

All lanes: HM13 Polyclonal Antibody at 6ug/ml+ Mouse liver tissue

Goat polyclonal to rabbit at 1/10000 dilution

Predicted band size: 41 kDa

Observed band size: 41 kDa



Immunofluorescent analysis of 293 cells using HM13 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)

*This product is for research purposes only. Not intended for use in diagnostic procedures.*

**EpiGentek Group Inc.** • USA Tel: 1-877-374-4368 • Int. Tel: 1-718-484-3990 • Fax: 1-718-484-3990 • [info@epigentek.com](mailto:info@epigentek.com) • [www.epigentek.com](http://www.epigentek.com)

2019-05-16-A59386