

EMP3 Polyclonal Antibody

(Catalog # A62470)

Background

Probably involved in cell proliferation and cell-cell interactions.

Description

EMP3 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

P54852

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Epithelial membrane protein 3 protein (25-65AA)

Storage

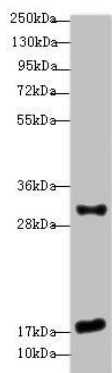
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Epithelial membrane protein 3, EMP-3, Hematopoietic neural membrane protein 1, HNMP-1, Protein YMP, EMP3, YMP

Application

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



Western Blot

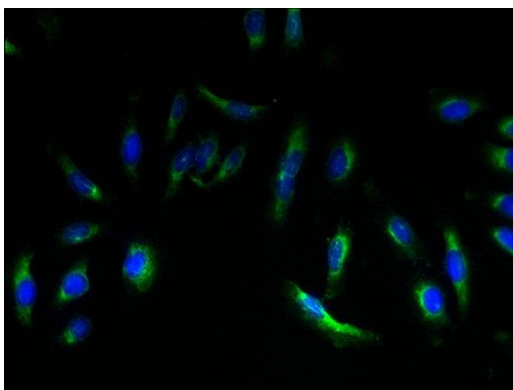
All lanes: EMP3 Polyclonal Antibody at 3 ug/ml+ Mouse heart tissue

Secondary

Goat polyclonal to rabbit IgG at 1/10000 dilution

Predicted band size: 18 kDa

Observed band size: 18, 31 kDa



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