
EZR Polyclonal Antibody

(Catalog # A71587)

Background

The cytoplasmic peripheral membrane protein encoded by this gene functions as a protein-tyrosine kinase substrate in microvilli. As a member of the ERM protein family, this protein serves as an intermediate between the plasma membrane and the actin cytoskeleton. This protein plays a key role in cell surface structure adhesion, migration and organization, and it has been implicated in various human cancers. A pseudogene located on chromosome 3 has been identified for this gene. Alternatively spliced variants have also been described for this gene.

Description

EZR 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

P15311

Purification

Affinity Purified

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 350-470 of human EZR (NP_003370.2)

Storage

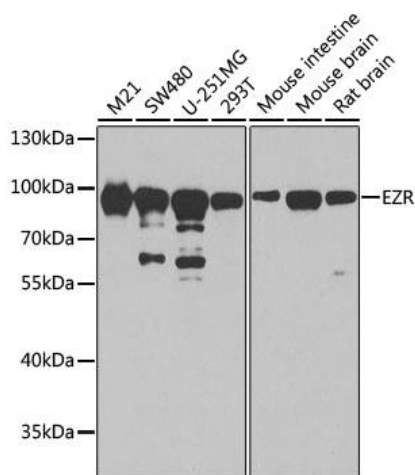
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

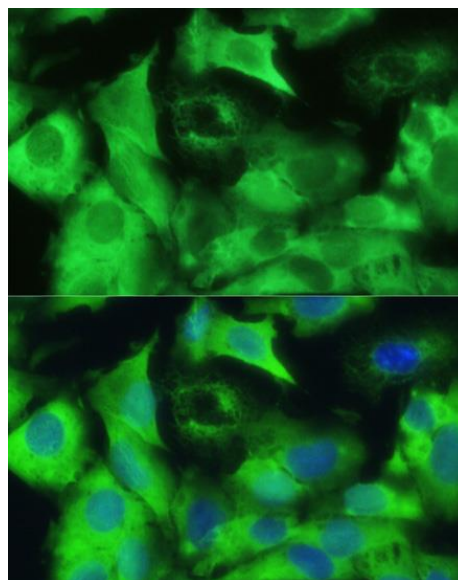
EZR; CVIL; CVL; HEL-S-105; VIL2; ezrin

Application

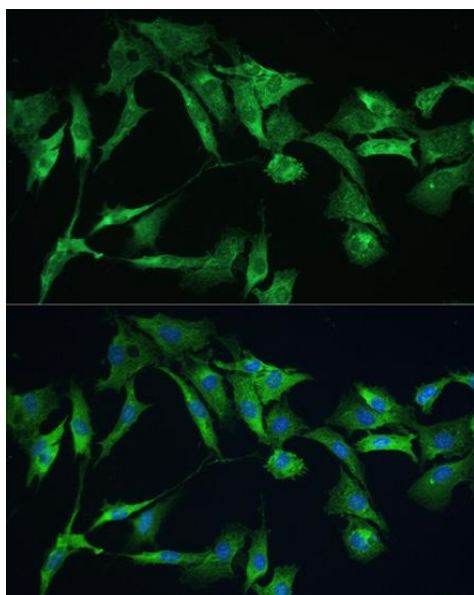
WB, IF, IP; Recommended dilution: WB 1:500 - 1:2000, IF/ICC 1:50 - 1:100, IP 1:50 - 1:100



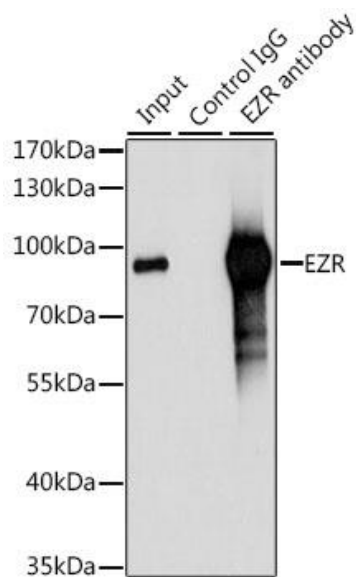
Western blot analysis of extracts of various cell lines, using EZR Polyclonal Antibody at 1:1000 dilution.
 Secondary Antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.
 Lysates/proteins: 25ug per lane.
 Blocking buffer: 3% nonfat dry milk in TBST.
 Exposure time: 30s.



Immunofluorescence analysis of U-2 OS cells using EZR Polyclonal Antibody at dilution of 1:100. Blue: DAPI for nuclear staining.



Immunofluorescence analysis of C6 cells using EZR Polyclonal Antibody at dilution of 1:100. Blue: DAPI for nuclear staining.



Immunoprecipitation analysis of 200ug extracts of 293T cells, using 3 ug EZR Polyclonal Antibody. Western blot was performed from the immunoprecipitate using EZR Polyclonal Antibody at a dilution of 1:1000.