

## TNFSF10 Monoclonal Antibody [2E1-1B9]

[Catalog No. A-6211-100]

### Description

Mouse monoclonal antibody raised against a full length recombinant TNFSF10, clone 2E1-1B9.

### Size

100 µg

### Purification

Protein A purified

### Immunogen

TNFSF10 (AAH32722, 1 a.a. ~ 282 a.a) full length recombinant protein with GST tag. MW of the GST tag alone is 26 KDa.

### Cross Reactivity

Human

### Isotype

IgG1 Kappa

### Formulation

In 1x PBS, pH 7.2

### Storage

Aliquot and store at -20°C or lower. Avoid repeated freezing and thawing. Multiple freeze/thaw cycles may result in decreased performance. Non-hazardous. No MSDS necessary.

### Handling Recommendations

For maximum recovery of the products, centrifuge the vial prior to opening the cap.

### Applications & Suggested Dilutions\*

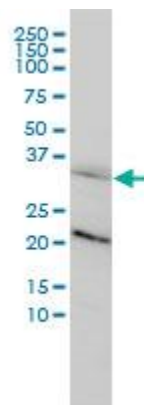
Western Blot: 0.5-1 µg/ml

ELISA: 0.25-1 µg/ml

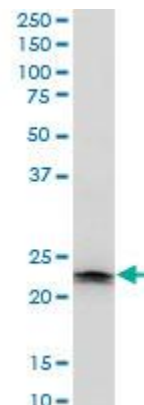
Immunohistochemistry (IHC): 2.5 µg/ml

\*The end user is responsible for determining optimal working dilutions.

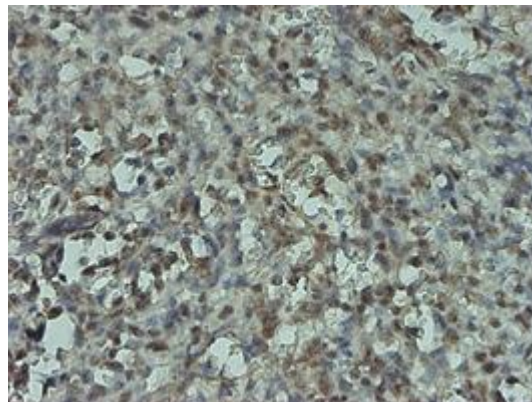
### Data



TNFSF10 Monoclonal Antibody, clone 2E1-1B9. Western Blot analysis of TNFSF10 expression in HeLa.



TNFSF10 Monoclonal Antibody, clone 2E1-1B9. Western Blot analysis of TNFSF10 expression in A-431.



Immunoperoxidase of monoclonal antibody to TNFSF10 on formalin-fixed paraffin-embedded human spleen. [Antibody Concentration: 2 µg/ml]

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