

## TRAF2 Monoclonal Antibody [4C11]

[Catalog No. A-6207-100]

### Description

Mouse monoclonal antibody raised against a full-length recombinant TRAF2, clone 4C11.

### Size

100 µg

### Purification

Protein A purified

### Immunogen

TRAF2 (AAH43492, 1 a.a. ~ 502 a.a) full-length recombinant protein with GST tag. MW of the GST tag alone is 26 KDa.

### Cross Reactivity

Human

### Isotype

IgG2a 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.

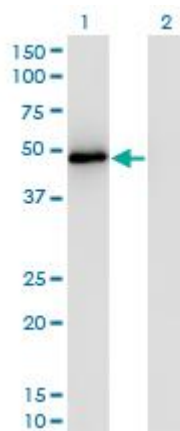
### Applications & Suggested Dilutions\*

Western Blot: 0.5-1 µg/ml

ELISA: 0.25-1 µg/ml

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

### Data



*Western Blot analysis of TRAF2 expression in transfected 293T cell line by TRAF2 Monoclonal Antibody, clone 4C11.*

*Lane 1: TRAF2 transfected lysate (55.9 KDa).*

*Lane 2: Non-transfected lysate.*

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