

## MAPK11 Monoclonal Antibody [1F9]

[Catalog No. A-7118-100]

### Description

Mouse monoclonal antibody raised against a partial recombinant MAPK11, clone 1F9.

### Size

100 µg

### Purification

Protein A purified

### Immunogen

MAPK11 (AAH27933, 255 a.a. ~ 364 a.a) partial 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.

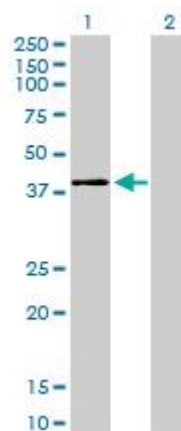
### Applications & Suggested Dilutions\*

Western Blot: 0.5-1 µg/ml

ELISA: 0.25-1 µg/ml

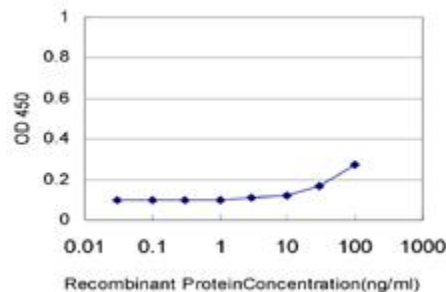
*\*The end user is responsible for determining optimal working dilutions.*

### Data



Western Blot analysis of MAPK11 expression in transfected 293T cell line by MAPK11 Monoclonal Antibody, clone 1F9.

Lane 1: MAPK11 transfected lysate (41.4 KDa).  
Lane 2: Non-transfected lysate.



Detection limit for recombinant GST tagged MAPK11 is approximately 10 ng/ml as a capture antibody.

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