

## MAP2K3 Monoclonal Antibody [1D10]

[Catalog No. A-7155-100]

### Description

Mouse monoclonal antibody raised against a partial recombinant MAP2K3, clone 1D10.

### Size

100 µg

### Purification

Protein A purified

### Immunogen

MAP2K3 (AAH32478, 1 a.a. ~ 100 a.a) partial 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\*

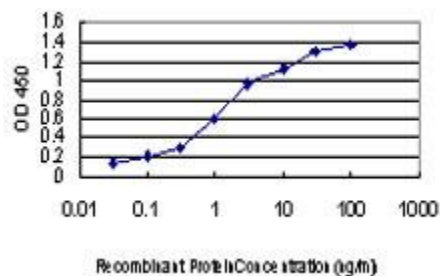
Western Blot: 0.5-1 µg/ml

ELISA: 0.25-1 µg/ml

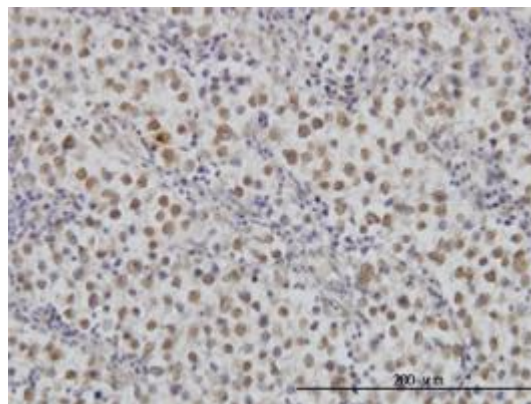
Immunohistochemistry (IHC): 2.5 µg/ml

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

### Data



*Detection limit for recombinant GST tagged MAP2K3 is approximately 0.03 ng/ml as a capture antibody.*



*Immunoperoxidase of monoclonal antibody to MAP2K3 on formalin-fixed paraffin-embedded in human dysgerminoma. [Antibody Concentration: 3 µg/ml]*

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