

## NEK2 Monoclonal Antibody [2F6]

[Catalog No. A-7139-100]

### Description

Mouse monoclonal antibody raised against a partial recombinant NEK2, clone 2F6.

### Size

100 µg

### Purification

Protein A purified

### Immunogen

NEK2 (AAH43502, 331 a.a. ~ 445 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\*

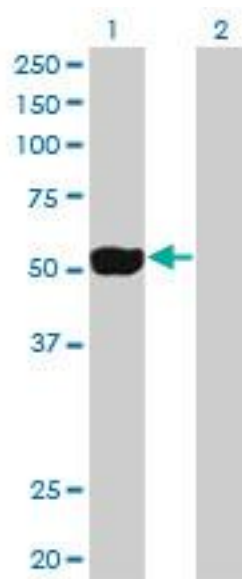
Western Blot: 0.5-1 µg/ml

ELISA: 0.25-1 µg/ml

Immunohistochemistry (IHC): 2.5 µg/ml

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

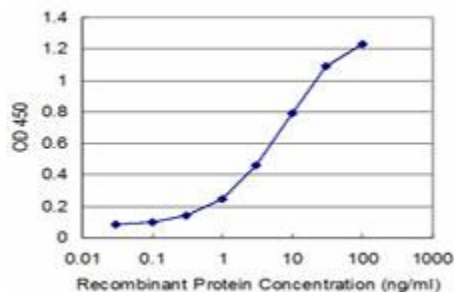
### Data



Western Blot analysis of NEK2 expression in transfected 293T cell line by NEK2 Monoclonal Antibody, clone 2F6.

Lane 1: NEK2 transfected lysate (51.8 KDa).

Lane 2: Non-transfected lysate.



Detection limit for recombinant GST tagged NEK2 is approximately 0.3 ng/ml as a capture antibody.

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