

## TLK1 Monoclonal Antibody [4B3]

[Catalog No. A-6124-100]

### Description

Mouse monoclonal antibody raised against a partial recombinant TLK1, clone 4B3.

### Size

100 µg

### Purification

Protein A purified

### Immunogen

TLK1 (AAH32657, 111 a.a. ~ 230 a.a) partial recombinant protein with GST tag. MW of the GST tag alone is 26 KDa.

### Cross Reactivity

Human

### Isotype

IgG3 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\*

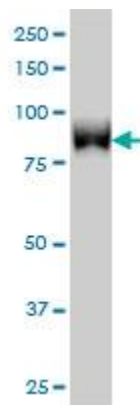
Western Blot: 0.5-1 µg/ml

ELISA: 0.25-1 µg/ml

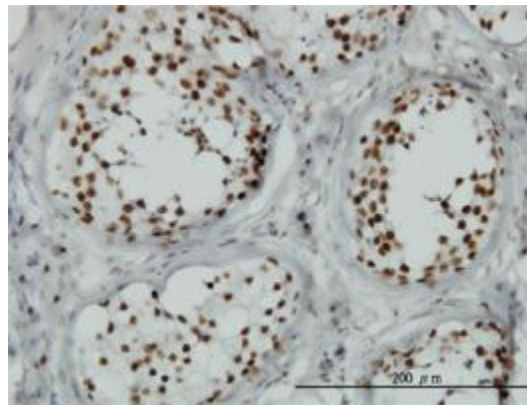
Immunohistochemistry (IHC): 2.5 µg/ml

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

### Data



TLK1 Monoclonal Antibody, clone 4B3. Western Blot analysis of TLK1 expression in HeLa NE.



Immunoperoxidase of monoclonal antibody to TLK1 on formalin-fixed paraffin-embedded human testis. [Antibody Concentration: 1 µg/ml]

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