

## CHEK2 Monoclonal Antibody [4B7]

[Catalog No. A-6123-100]

### Description

Mouse monoclonal antibody raised against a full length recombinant CHEK2, clone 4B7.

### Size

100 µg

### Purification

Protein A purified

### Immunogen

CHEK2 (AAH04207, 1 a.a. ~ 544 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

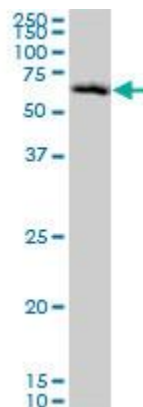
ELISA: 0.25-1 µg/ml

Immunohistochemistry (IHC): 2.5 µg/ml

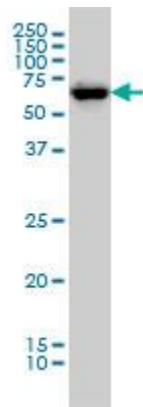
Immunofluorescence: 5 µg/ml

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

### Data



*CHEK2 Monoclonal Antibody, clone 4B7. Western Blot analysis of CHEK2 expression in HeLa.*



*CHEK2 Monoclonal Antibody, clone 4B7. Western Blot analysis of CHEK2 expression in HeLa NE.*

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