

## UBE2M Monoclonal Antibody [3C4]

[Catalog No. A-8114-100]

### Description

Mouse monoclonal antibody raised against a full length recombinant UBE2M, clone 3C4.

### Size

100 µg

### Purification

Protein A purified

### Immunogen

UBE2M (AAH58924, 1 a.a. ~ 184 a.a) full length 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\*

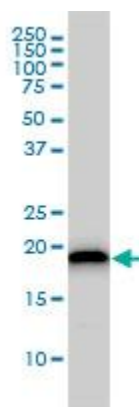
Western Blot: 0.5-1 µg/ml

ELISA: 0.25-1 µg/ml

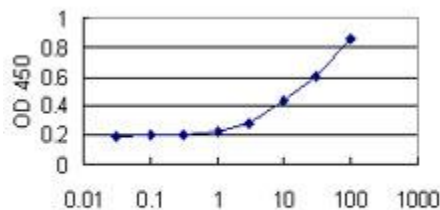
Immunofluorescence: 5 µg/ml

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

### Data

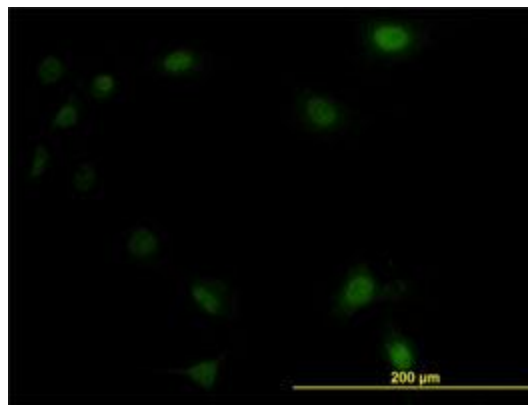


UBE2M Monoclonal Antibody, clone 3C4. Western Blot analysis of UBE2M expression in Jurkat.



Recombinant Protein Concentration (ng/ml)

Detection limit for recombinant GST tagged UBE2M is approximately 0.1 ng/ml as a capture antibody.



Immunofluorescence of monoclonal antibody to UBE2M on HeLa cell. [Antibody Concentration: 10 µg/ml]

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