

## UHRF2 Monoclonal Antibody [3A11]

[Catalog No. A-8139-100]

### Description

Mouse monoclonal antibody raised against a partial recombinant UHRF2, clone 3A11.

### Size

100 µg

### Purification

Protein A purified

### Immunogen

UHRF2 (NP\_690856, 81 a.a. ~ 181 a.a) partial recombinant protein with GST tag. MW of the GST tag alone is 26 KDa.

### Cross Reactivity

Human

### Isotype

IgG2b 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.

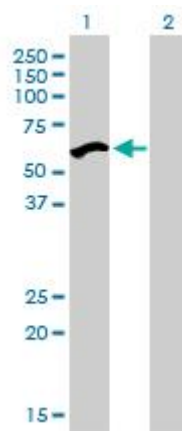
### Applications & Suggested Dilutions\*

Western Blot: 0.5-1 µg/ml

ELISA: 0.25-1 µg/ml

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

### Data



*Western Blot analysis of UHRF2 expression in transfected 293T cell line by UHRF2 Monoclonal Antibody, clone 3A11.*

*Lane 1: UHRF2 transfected lysate (56.1 KDa).*

*Lane 2: Non-transfected lysate.*

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