

## SENP8 Monoclonal Antibody [2E1]

[Catalog No. A-8149-100]

### Description

Mouse monoclonal antibody raised against a partial recombinant SENP8, clone 2E1.

### Size

100 µg

### Purification

Protein A purified

### Immunogen

SENP8 (NP\_660205, 113 a.a. ~ 213 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.

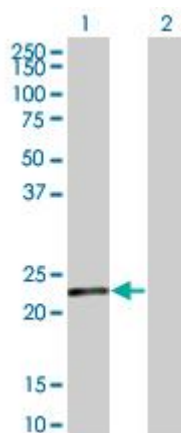
### Applications & Suggested Dilutions\*

Western Blot: 0.5-1 µg/ml

ELISA: 0.25-1 µg/ml

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

### Data



Western Blot analysis of SENP8 expression in transfected 293T cell line by SENP8 Monoclonal Antibody, clone 2E1.

Lane 1: SENP8 transfected lysate (24.1 KDa).  
Lane 2: Non-transfected lysate.

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