

SENP6 Monoclonal Antibody [4B7]

[Catalog No. A-8147-100]

Description

Mouse monoclonal antibody raised against a partial recombinant SENP6, clone 4B7.

Size

100 µg

Purification

Protein A purified

Immunogen

SENP6 (NP_056386, 1 a.a. ~ 111 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*

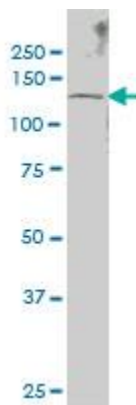
Western Blot: 0.5-1 µg/ml

ELISA: 0.25-1 µg/ml

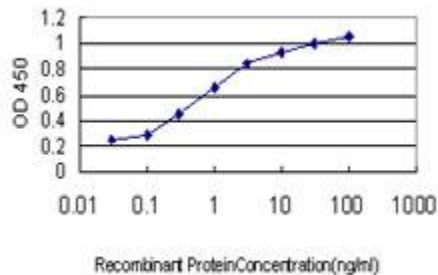
Immunohistochemistry (IHC): 5 µg/ml

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

Data



SENP6 Monoclonal Antibody, clone 4B7. Western Blot analysis of SENP6 expression in IMR-32.



Detection limit for recombinant GST tagged SENP6 is approximately 0.03 ng/ml as a capture antibody.

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