

FADD Monoclonal Antibody [3A12]

[Catalog No. A-6221-100]

Description

Mouse monoclonal antibody raised against a partial recombinant FADD, clone 3A12.

Size

100 µg

Purification

Protein A purified

Immunogen

FADD (AAH00334, 109 a.a. ~ 208 a.a) partial recombinant protein with GST tag. MW of the GST tag alone is 26 KDa.

Cross Reactivity

Human

Isotype

IgG1 Lambda

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.

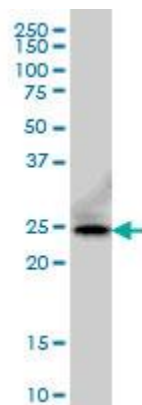
Applications & Suggested Dilutions*

Western Blot: 0.5-1 µg/ml

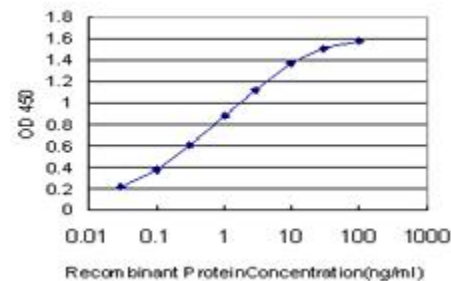
ELISA: 0.25-1 µg/ml

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

Data



FADD Monoclonal Antibody, clone 3A12. Western Blot analysis of FADD expression in A-431.



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

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