

MAPK13 Monoclonal Antibody [2D8]

[Catalog No. A-7120-100]

Description

Mouse monoclonal antibody raised against a partial recombinant MAPK13, clone 2D8.

Size

100 µg

Purification

Protein A purified

Immunogen

MAPK13 (AAH00433, 251 a.a. ~ 365 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

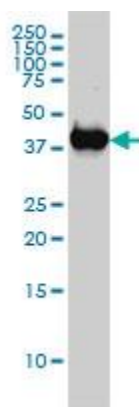
ELISA: 0.25-1 µg/ml

Immunohistochemistry (IHC): 2.5 µg/ml

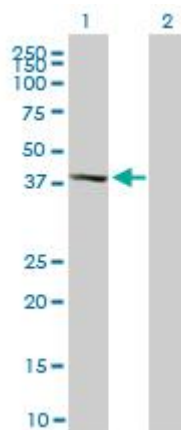
Immunofluorescence: 5 µg/ml

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

Data



MAPK13 Monoclonal Antibody, clone 2D8. Western Blot analysis of MAPK13 expression in MCF-7.



Western Blot analysis of MAPK13 expression in transfected 293T cell line by MAPK13 Monoclonal Antibody, clone 2D8.

Lane 1: MAPK13 transfected lysate (42.1 KDa).
Lane 2: Non-transfected lysate.

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