

[pSer218/222]MEK1 and [pSer222/226]MEK2 monoclonal antibody (7E10)

MEK is a dual specificity kinase capable of phosphorylating both tyrosine and threonine residues. The MEK family, which is also known as MAP kinase kinase, phosphorylates MAP kinases on the conserved T-X-Y motif; phosphorylation of MAPK by MEK results in an increase in MAPK activity. MEK is involved in a diverse array of cellular processes such as stress-activated response, apoptosis, cytokine-induced cell proliferation, and DNA recombination during meiosis.

This antibody is covered by our [Worry-Free Guarantee](#).

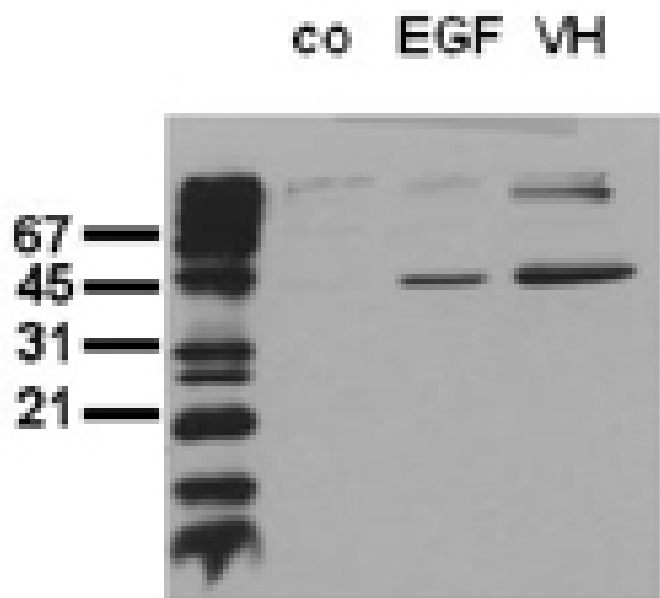
Ordering Information

[Order Online »](#)

ADI-905-652-100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of untreated control (co), EGF stimulated (EGF), or pervanadate treated (VH) A549 cell lysates, probed with [pS218/222]MEK1 [pS222/226]MEK2, mAb (7E10) (0.5 μ g/ml).

Handling & Storage

Use/Stability Stable at -80°C up to 1 year, at 4°C up to 3 months.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name MKK1/MKK2, MAP kinase kinase 1/2

Application ELISA, WB

Application Notes Detects a band of ~45kDa by Western blot.

Clone 7E10

Formulation Lyophilized from 1ml of 2x PBS containing 0.09% sodium azide, PEG, and sucrose.

Host Mouse

Immunogen Synthetic peptide corresponding to a portion of MEK1 phosphorylated at Ser218/222 and MEK2 phosphorylated at Ser222/226 , conjugated to KLH.

Isotype IgG1

Purity Detail Thiophilic adsorption and size exclusion chromatography purified.

Recommendation Dilutions/Conditions ELISA (0.1µg/ml)Western Blot (0.5µg/ml, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Reconstitution Reconstitute with 1ml water (15 minutes at room temperature).

Species Reactivity Dog, Human, Mouse

UniProt ID

Q02750 (MEK1), P36507 (MEK2)

Worry-free Guarantee

This antibody is covered by our [Worry-Free Guarantee](#)

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)