# [pSer<sup>218/222</sup>]MEK1 and [pSer<sup>222/226</sup> ]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**

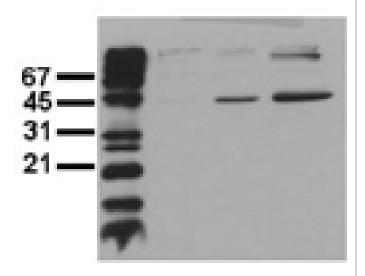
Order Online »

**ADI-905-652-100** 100μg

Manuals, SDS & CofA

**View Online »** 

## co EGF VH



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 µg/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 Ser218/222and 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

**Worry-free Guarantee** 

This antibody is covered by our Worry-Free Guarantee

.

Last modified: May 29, 2024



 ENZO LIFE SCIENCES,
 European Sales Office

 INC.
 ENZO LIFE SCIENCES

 Phone: 800.942.0430
 (ELS) AG

 info Phone: +41 61 926 8989

 usa@enzolifesciences.com
 info

eu@enzolifesciences.com

Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany Phone: +49 7621 5500 526 infode@enzolifesciences.com UK & Ireland Phone (UK customers): 0845 601 1488 Phone: +44 1392 825900 infouk@enzolifesciences.com