

# ERK1/2 monoclonal antibody (12D4)

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

Citations: 1

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-804-061-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

# Handling & Storage

Use/Stability	Stable at -80°C up to 1 year, at 4°C up to 3 months.
Handling	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and freeze in liquid nitrogen.
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

# Product Details

Alternative Name	MAPK2/3, MAP kinase 2/3
Application	ELISA, ICC, IHC, IP, WB
Clone	12D4
Crossreactivity	Does not cross-react with activated JNK or p38.
Formulation	Lyophilized from 1ml of 2x PBS containing 0.09% sodium azide, PEG, and sucrose.
Host	Mouse
Immunogen	Peptide comprising the pThr-Glu-pTyr motif in activated ERK1 and ERK2 conjugated to KLH.
Isotype	IgG1
Positive Control	Included. (Prod. No. ALX-840-025)
Purity Detail	Thiophilic adsorption and size exclusion chromatography purified.
Recommendation Dilutions/Conditions	Immunocytochemistry (1-10µg/ml)Immunoprecipitation (1-10µg per 106 pervanadate-treated A431 or HepG2 cells)Western Blot (0.5µg/ml for HRPO/ECL detection; recommended blocking buffer CPPT: 10mM TRIS-HCl, pH 7.4, 0.5% (w/v) casein, 1% (w/v) PEG 4,000, 1% (w/v) polyvinylpyrrolidone, 0.1% (v/v) Tween 20, 150mM sodium chloride)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).

<b>Species Reactivity</b>	Dog, Human, Mouse, Rat
<b>Specificity</b>	Recognizes the pThr-Glu-pTyr motif of activated ERK1 and ERK2. Requires phosphorylation both at the threonine and the tyrosine site and does not cross-react with the non-phosphorylated form of the protein.
<b>Technical Info / Product Notes</b>	Replacement for <b>ADI-905-651</b>
<b>UniProt ID</b>	P27361 (Erk1), P28482 (Erk2)
<b>Worry-free Guarantee</b>	This antibody is covered by our <a href="#">Worry-Free Guarantee</a> .



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)