

Erk1/2 polyclonal antibody

The Extracellular regulated kinases, Erk-1 and Erk-2, both play a critical role in cellular proliferation. These extracellular kinases, are known to reside downstream of the Mitogen activated protein kinase/Extracellular protein kinases referred to as MEK. Both MEK-1 and MEK-2 phosphorylate and activate Erk on the conserved TEY motif within the activation loop of the kinase.

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

Citations: 14

[View Online »](#)

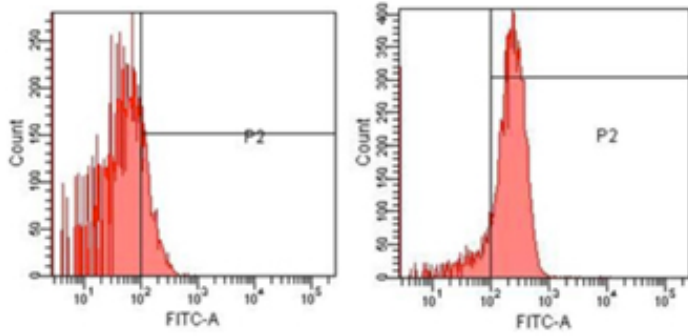
Ordering Information

[Order Online »](#)

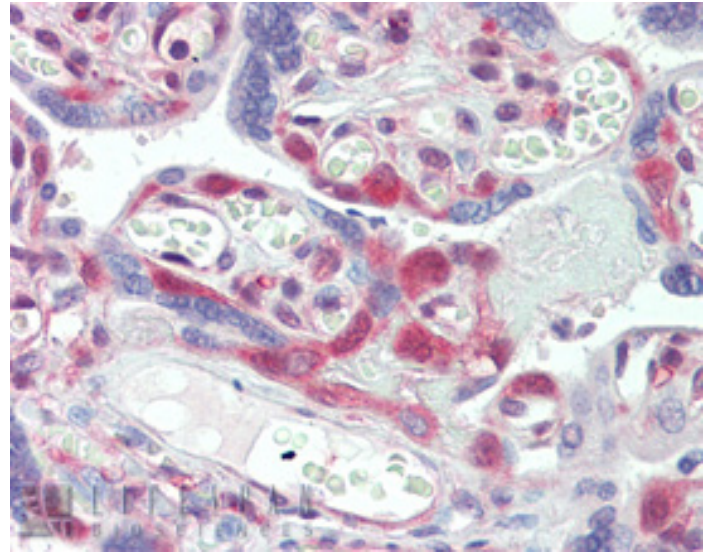
ADI-KAP-MA001-D	50µg
ADI-KAP-MA001-F	200µg

Manuals, SDS & CofA

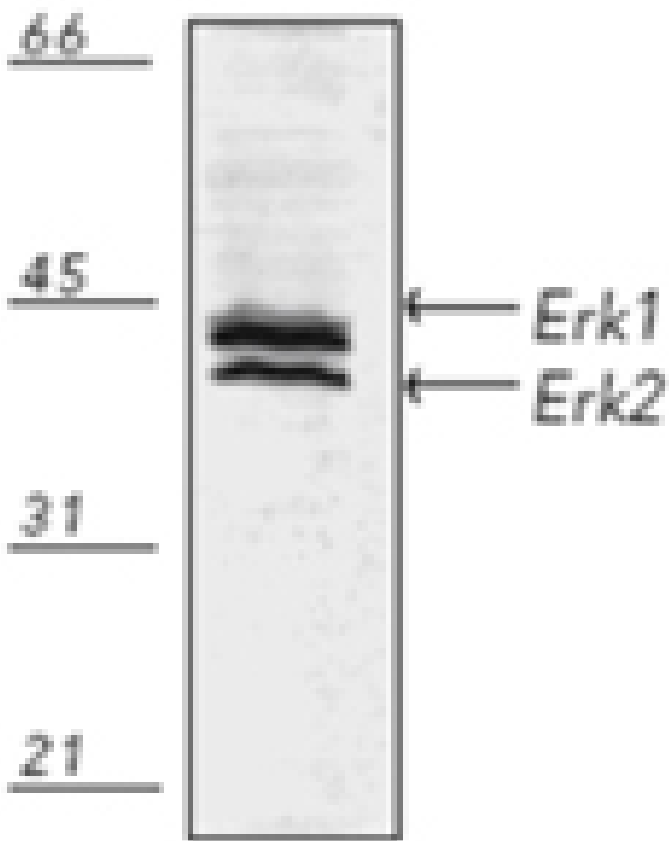
[View Online »](#)



Flow cytometry analysis of U251 cells using Erk1/2 pAb (right) and control rabbit IgG (left).



Immunohistochemistry analysis: Human placenta tissue stained with Erk1/2, pAb at 10µg/ml.



Western blot analysis of human A431 fibroblast cell lysate, probed with Erk1/2 pAb.

Handling & Storage

Handling Avoid freeze/thaw cycles.

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 Flow Cytometry, IHC (PS), IP, WB

Application Notes Detects a bands of ~42kDa (Erk2) and ~44kDa (Erk1) by Western blot.

Formulation Liquid. In PBS containing 50% glycerol and 0.09% sodium azide.

GenBank ID M61177

Host Rabbit

Immunogen Synthetic peptide corresponding to a portion of rat Erk1. The sequence is completely conserved in hamster.

Purity Detail Protein A affinity purified.

Recommendation Flow Cytometry (10µg/ml)Immunohistochemistry (paraffin sections, 1:50)Western Blot
Dilutions/Conditions (1:1,000, colorimetric)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Source Purified from rabbit serum.

Species Reactivity Bovine, Chicken, Drosophila, Human, Mouse, Rat, Sheep, Xenopus

UniProt ID P21708

Worry-free Guarantee

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



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)