

[pThr185/pTyr187]ERK1/2 polyclonal antibody

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

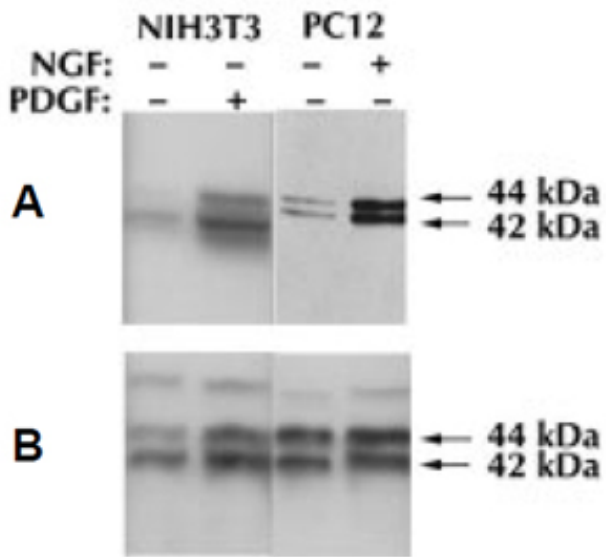
Ordering Information

[Order Online »](#)

BML-SA275-0020	20µl
----------------	------

Manuals, SDS & CofA

[View Online »](#)



A. ERK1 & 2 (pTpY 185/187). B. ERK1 & 2.
 PC12 cells were treated with NGF and NIH3T3 cells were treated with PDGF and cell extracts were immunoblotted with either [pThr¹⁸⁵/pTyr¹⁸⁷]ERK1/2, pAb (Prod. No. BML-SA275) or ERK1/2, pAb (Prod. No. BML-SA278).

Handling & Storage

Long Term Storage -20°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	MAPK2/3, MAP kinase 2/3
Application	IHC, WB
Formulation	Liquid. In Dulbecco's PBS, pH 7.3, containing 50% glycerol, 1.0 mg/ml BSA, and 0.05% sodium azide.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to a portion of human ERK1&2 phosphorylated at Thr202/185 and Tyr204/187. This region is conserved among many species including mouse, rat, bovine, frog and fly.
Purity Detail	Purified by sequential epitope-specific chromatography. The antibody has been negatively preadsorbed using a non-phosphopeptide corresponding to the phosphopeptide used as the immunogen. The final product was generated by affinity chromatography.
Species Reactivity	Chicken, Human, Mouse, Rat
Specificity	Recognizes ERK1/2 phosphorylated at pThr ¹⁸⁵ /Tyr ¹⁸⁷ .
UniProt ID	P27361 (Erk1), P28482 (Erk2)
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

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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