

ERK1/ERK2 (172-190)

MEK substrate

This sequence, ADPDHDHTGFLTGYVATRRR, is based on Erk1/Erk2 residues 172-190, containing a MAP kinase kinase phosphorylation site (underlined). N-terminal Arg is added for phosphocellulose binding.

Ordering Information

[Order Online »](#)

BML-P124-0001	1mg
----------------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

MW	2543.8
Purity	≥97%
Sequence	Ala-Asp-Pro-Asp-His-Asp-His-Thr-Gly-Phe-Leu-Thr-Glu-Tyr-Val-Ala-Thr-Arg-Trp-Arg-Arg

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