

[pThr183, pTyr185] MAP kinase (177-189)

VHR substrate

Diphosphopeptide with sequence, DHTGFLpTEpYVATR, corresponding to MAP kinase residues 177-189 inclusive of putative activation sites. Peptide can be used as a VHR phosphatase substrate (BML-SE333). Phosphate release may be quantified using BML-AK111 malachite green colorimetric reagent. Phosphate release is much more rapid for the phosphotyrosine residue compared to the phosphothreonine residue, both *in vitro* and *in vivo*.

Ordering Information

[Order Online »](#)

BML-P164-0001	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

Appearance	White powder.
MW	1669.6
Purity	≥95% (HPLC)
Sequence	Asp-His-Thr-Gly-Phe-Leu-pThr-Glu-pTyr-Val-Ala-Thr-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