

# [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*.

**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](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)