

R18 peptide

14-3-3 inhibitor

The peptide sequence, PHCVPRDLSWLDLEANMCLP, named R18, was identified as an unphosphorylated peptide that binds to 14-3-3 proteins with high affinity ($K_d=80$ nM). It has been shown to bind to the same amphipathic groove on the surface of 14-3-3 as phosphorylated peptides. R18 can inhibit the formation of the 14-3-3-Raf-1 complex, reduce Raf-1 dependent transcriptional activity in cells and block the activation of ExoS by 14-3-3. It binds to all the 14-3-3 isoforms with equal affinities. This peptide is useful for competitive inhibition experiments. A biotinylated form is available (BML-P225), allowing detection using streptavidin.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-P214-0001	1mg
---------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Solid.
Formulation	Lyophilized. Trifluoroacetate salt.
Identity	Determined by MS.
MW	2309.6
Purity	≥95% (HPLC)
Sequence	Pro-His-Cys-Val-Pro-Arg-Asp-Leu-Ser-Trp-Leu-Asp-Leu-Glu-Ala-Asn-Met-Cys-Leu-Pro-OH
Solubility	Soluble in Water.
Source	Synthetic.

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-fr@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