

# R18 peptide (biotinylated)

## 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 and the N-terminal biotin allows it to be used in assay with streptavidin bound membranes.

### Ordering Information

[Order Online »](#)

BML-P225-0001	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** Lyophilized solid.

**MW** 2536.7

**Purity** ≥95%

**Sequence** Biotin-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.



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)