

Crosstide (biotinylated)

Akt substrate

This peptide sequence, biotin-GRPRTSSFAEG, corresponds to the sequence of GSK3 surrounding the serine phosphorylated by MAP kinase activated protein kinase (MAPKAP kinase-1) and p70 ribosomal S6 kinase. Acts as a substrate for Akt (PKB), $K_m=4 \mu\text{M}$. The N-terminal biotin allows this peptide to be used in kinase assays with streptavidin-bound membranes.

Ordering Information

[Order Online »](#)

BML-P191-0001	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	1390.4
Purity	≥95%
Sequence	Biotin-Gly-Arg-Pro-Arg-Thr-Ser-Ser-Phe-Ala-Glu-Gly
Solubility	Soluble in Water.

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