

Crosstide

Akt substrate

This peptide sequence, 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), $K_m=4\mu\text{M}$. Useful in phosphocellulose kinase assays.

Ordering Information

[Order Online »](#)

BML-P149-0001	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

CAS 171783-05-4

Formula $C_{48}H_{77}N_{17}O_{17}$

MW 1165.2

Sequence Gly-Arg-Pro-Arg-Thr-Ser-Ser-Phe-Ala-Glu-Gly

Technical Info / Product Notes Replacement for **ADI-SPK-108**

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-
fr@enzolifesciences.com](mailto:info-fr@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)