

GSK3 substrate

This peptide sequence, RRRAAEELDSRAGpSPQL, corresponds to the sequence of the ϵ subunit of eIF2B surrounding Ser⁵⁴⁰, a target of GSK-3 (glycogen synthase kinase 3) phosphorylation, with the addition of a phosphoserine in the 4 position. This phosphoserine allows for more efficient phosphorylation of the second serine in the peptide (K_m = 49 μ M for GSK-3 β , 83 μ M for GSK-3 α). Useful in phosphocellulose GSK-3 α and - β assays.

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-P151-0001	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

Alternative Name Glycogen synthase kinase 3 substrate

MW 1992.0

Sequence H-Arg-Arg-Arg-Ala-Ala-Glu-Glu-Leu-Asp-**Ser**-Arg-Ala-Gly-**pSer**-Pro-Gln-Leu-OH

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)