

Ribosomal protein S6 (rat) (229-239)

S6 kinase substrate

This sequence, AKRRRLSSLRA, derived from rat hepatic 40S ribosomal protein S6 (229-239). Substrate for Rho-kinase II (ROCK II), $K_m=0.9\text{ }\mu\text{M}$ and for p90 rsk (S6K) in cell extracts ($K_m=0.18\text{ mM}$) using phosphocellulose assay. Peptide phosphorylation by PKS and PKC can be eliminated by the inclusion of PKI and EGTA, respectively, in the assay.

Ordering Information

[Order Online »](#)

BML-P144-0001	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

MW 1313.9

Sequence Ala-Lys-Arg-Arg-Arg-Leu-Ser-Ser-Leu-Arg-Ala

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)