

SOS-1 (human) (1149-1158)

PI3 kinase inhibitor

SH3 binding sequence of guanine nucleotide exchange factor hSOS which is a ligand for the adaptor protein Grb2 SH3 domain. Blocks SOS/Grb2 interaction and thereby prevents Ras activation via receptor protein tyrosine kinases. $K_m = 5.6 \mu M$ for N-terminal SH3 of Grb2.

Ordering Information

[Order Online »](#)

BML-P305-0001	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 SOS-1 (human) N10 SH3 binding domain, Ras inhibitory peptide

Appearance Lyophilized solid.

MW 1169.7

Purity ≥98%

Sequence Val-Pro-Pro-Pro-Val-Pro-Pro-Arg-Arg-Arg

Solubility Soluble in Water.

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)