

pp60^{c-src} C-terminal phosphoregulatory peptide

Src kinase inhibitor

This phosphopeptide is the C-terminal negative regulatory domain of c-Src (521-533) which includes the negative regulatory site (Tyr⁵²⁷). This peptide, when tyrosine-phosphorylated, binds to the internal SH2 domain.

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-P301-0001	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 [pTyr527] c-Src (521-533)

Appearance Lyophilized solid.

MW 1585.7

Purity ≥98%

Sequence Ac-Thr-Ser-Thr-Glu-Pro-Gln-**pTyr**-Gln-Pro-Gly-Glu-Asn-Leu

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)