

Peptide A

Src kinase inhibitor

Sequence is pp60^{v-src} (137-157), which is located in the SH3/SH2 domain region. Non-competitive inhibitor of Src and EGFR. Selectively inhibits protein tyrosine phosphorylation in cell extracts. Peptide is not phosphorylated by Src. $IC_{50}=7.5\text{ }\mu\text{M}$ for v-Src. Inhibitory target seems to be a catalytic negative regulatory domain of the tyrosine kinase.

Ordering Information

[Order Online »](#)

BML-P306-0500	0.5mg
---------------	-------

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 pp60v-src (137-157)

MW 2482.8

Purity ≥97%

Sequence Val-Ala-Pro-Ser-Asp-Ser-Ile-Gln-Ala-Glu-Glu-Trp-Tyr-Phe-Gly-Lys-Ile-Thr-Arg-Arg-Glu



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)