

pp60^{v-src}

autophosphorylation site (phosphorylated)

PTP substrate

This sequence, RRLIEDNEpYTARG, is the v-Src (412-422) autophosphorylation site and this peptide contains a phosphorylated tyrosine. Useful in a nonradioactive tyrosine phosphatase assay and as a competitive tyrosine phosphatase inhibitor.

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 1671.9

Sequence Arg-Arg-Leu-Ile-Glu-Asp-Asn-Glu-**pTyr**-Thr-Ala-Arg-Gly

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com