

[pTyr1146/1150]

Insulin receptor (1142-1153)

Insulin receptor substrate

This sequence, TRDlpYETDpYYRK, is from the insulin receptor β -subunit cytoplasmic domain, inclusive of regulatory autophosphorylation sites Tyr¹¹⁴⁶, Tyr¹¹⁵⁰ and Tyr¹¹⁵¹. The unphosphorylated version of this peptide has been used as an insulin receptor kinase substrate. The phosphorylated peptide has been used as a protein tyrosine phosphatase substrate.

Ordering Information

[Order Online »](#)

BML-P321-0500	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

MW	1782.6
Sequence	Thr-Arg-Asp-Ile- pTyr -Glu-Thr-Asp- pTyr -Tyr-Arg-Lys

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)