

IRS-1 (Tyr608) peptide

Insulin receptor substrate

This sequence, KKHTDDGYMPMSPGVA, is from insulin receptor substrate-1 (IRS-1) inclusive of Tyr⁶⁰⁸ (mouse) -Tyr⁶¹² (human). It contains the insulin receptor tyrosine kinase substrate motif YMXM (Tyr-Met-X-Met). This peptide has been used as a substrate for purified insulin receptor ($K_m = 90 \mu M$) and other tyrosine kinases in phosphocellulose binding assays. The tyrosine phosphorylated version of this peptide binds to phosphatidylinositol 3-kinase (PI 3-kinase) SH2 domain and activates the enzyme.

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

| | |
|---------------|-----|
| BML-P320-0001 | 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 Insulin receptor substrate (Tyr608) peptide

Appearance White to off-white powder.

Formulation Lyophilized.

MW 1732.8

Purity ≥95% (HPLC)

Sequence Lys-Lys-His-Thr-Asp-Asp-Gly-Tyr-Met-Pro-Met-Ser-Pro-Gly-Val-Ala

Solubility Soluble in water.

Source Synthetic.

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)