

[pTyr992] EGF receptor (988-998)

PTP1B substrate

This sequence, DADEpYLIPQQG, is from an autophosphorylation site (Tyr⁹⁹²) of epidermal growth factor receptor (EGFR). An excellent substrate for mammalian PTP1B with $K_m = 3.9 \mu\text{M}$) and *Yersinia* PTPase. May be used in non-radioactive phosphatase assays incorporating the BIOMOL GREENTM reagent.

Ordering Information

[Order Online »](#)

BML-P323-0001	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 PTP1B substrate, [pTyr992] EGF receptor (988-998)]

Appearance White to off-white powder.

Formulation Lyophilized.

Identity Determined by MS.

MW 1330.3

Purity ≥95% (HPLC)

Sequence Asp-Ala-Asp-Glu-**pTyr**-Leu-Ile-Pro-Gln-Gln-Gly

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-
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)