

PTP1B (human), (recombinant)

PTP1B is the ubiquitous, prototypical nontransmembrane PTP originally identified in human placenta.

Citations: 36

[View Online »](#)

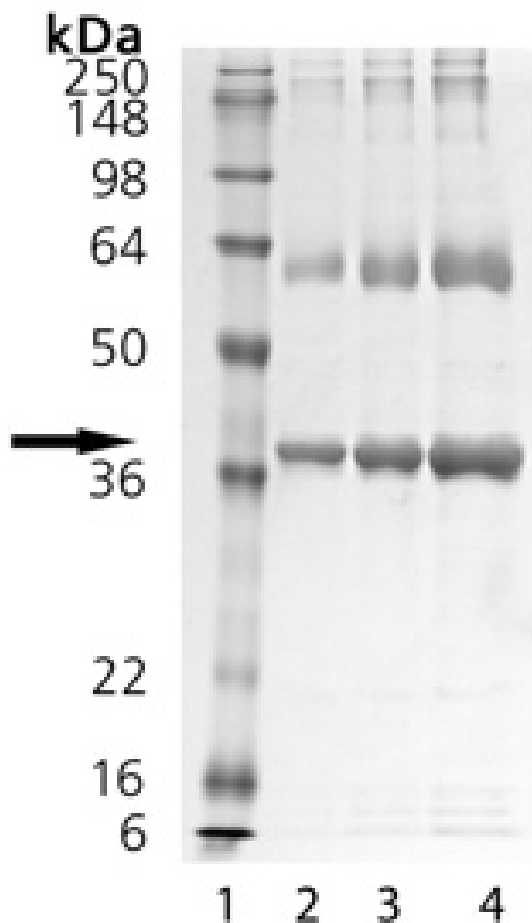
Ordering Information

[Order Online »](#)

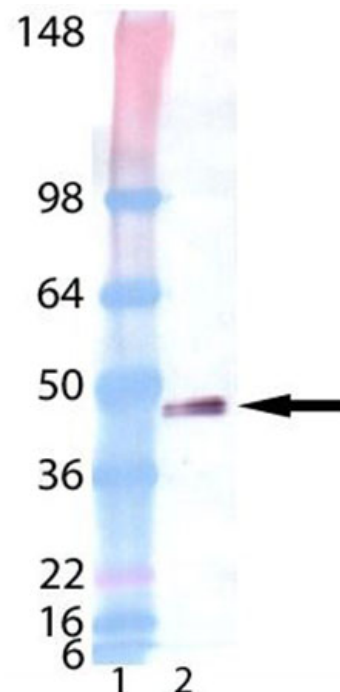
BML-SE332-0050	50µg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE Analysis: Lane 1: MW Marker, Lane 2: 1 µg, Lane 3: 2 µg, Lane 4: 5 µg PTP1B. The higher molecular weight bands are BSA included for stability in storage buffer.



Western Blot analysis of PTP1B (human), (recombinant) (Prod. No. BML-SE332): Lane 1: MW Marker, Lane 2: 100ng of PTP1B (human), (recombinant) probed with Anti-6X His tag® antibody.

Handling & Storage

Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -80°C.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Protein tyrosine phosphatase 1B
Application Notes	Useful for the study of tyrosine phosphatase kinetics, substrate specificity and for screening inhibitors.
Formulation	Liquid. In 50mM HEPES, pH7.2, containing 1mM EDTA, 5mM DTT, and 0.05% NP-40.
MW	~37.4kDa
Purity	≥95% (SDS-PAGE)
Purity Detail	Purified by multi-step chromatography.
Source	Produced in <i>E. coli</i> . Human PTP1B (protein tyrosine phosphatase 1B) (aa 1-322).
Specific Activity	≥20 μmol/min/mg at 30°C using 0.15mM of phosphopeptide substrate based on EGFR residues 988-998 (Prod. No. BML-P323) in 50mM HEPES, pH 7.2, 1mM EDTA, 1mM DTT, 0.05% NP-40, and 0.2μg/μl BSA.
UniProt ID	P18031

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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@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)