

T cell protein tyrosine phosphatase (human), (recombinant)

Citations: 3

[View Online »](#)

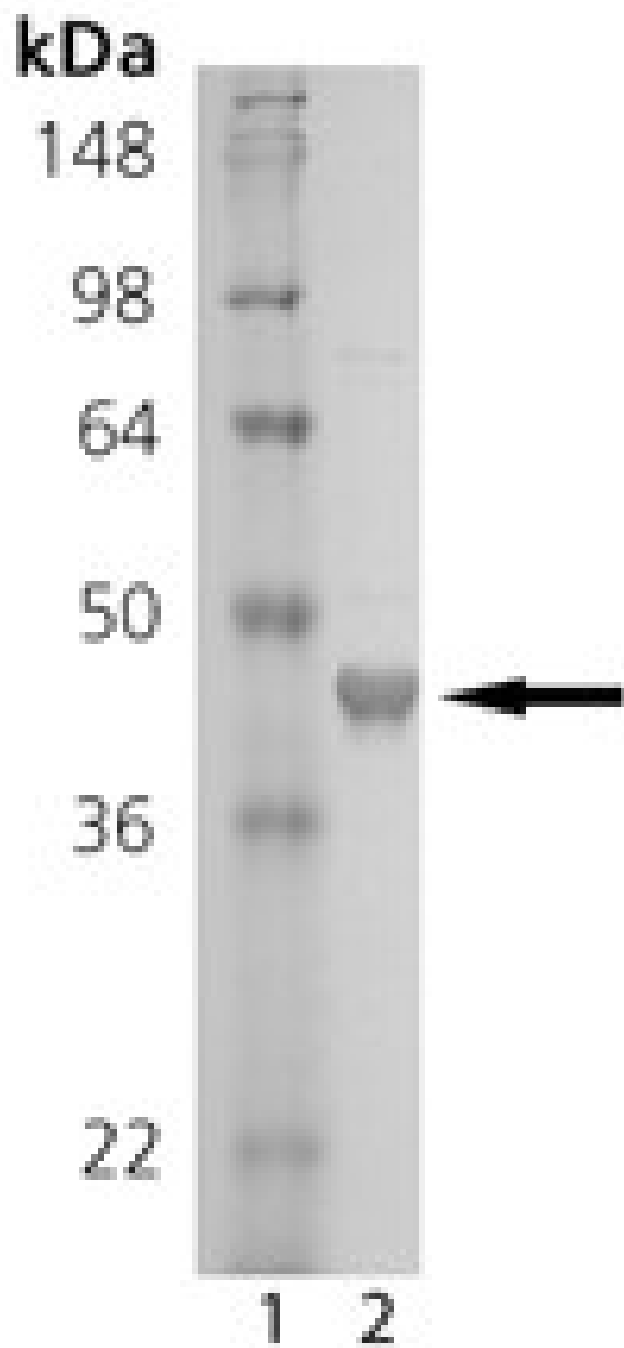
Ordering Information

[Order Online »](#)

BML-SE114-0200	200U
----------------	------

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE Analysis: Lane 1: MWM; Lane 2: 0.5 μ g of purified T-cell protein tyrosine phosphatase.

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	TCPTP, T cell PTP
Formulation	Liquid. In 50mM Tris-HCl, pH 7.0 containing 100 mM NaCl, 2 mM EDTA, 5 mM DTT, 0.01% Brij35 and 30% glycerol.
MW	38 kDa
Purity	≥95% (SDS-PAGE)
Purity Detail	Contains no detectable protease, DNase, RNase, or Ser/Thr phosphatase activity.
Source	Produced in <i>E. coli</i> . T-cell phosphatase with a C-terminal 11 kDa deletion (TCΔC11) (aa 1-317).
Specific Activity	One unit will release 1 nmol p-nitrophenol per minute at 30°C from 50mM p-nitrophenyl phosphate in 25mM TRIS/HCl, 50mM NaCl, 2mM EDTA, 5mM DTT, 0.01% Brij35, 1mg/ml BSA, pH 7.0.
UniProt ID	P17706

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)