

PTX3 (human), (recombinant)

Citations: 11

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-201-124-C010	10µg
------------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Pentraxin 3, TSG14, TNF stimulated gene-14

Application Notes ELISA and Western blot standard together with MAb to PTX3 (Prod. No. ALX-804-464 (capture antibody)) and PAb to PTX3-Biotin (Prod. No. ALX-210-365B (detection antibody)).

Concentration 0.1mg/ml.

Formulation Liquid. In PBS containing 2% BSA and 0.1% sodium azide.

MW ~42kDa. Due to glycosylation, PTX3 may migrate between 40-45kDa.

Source Produced in CHO cells. Human full-length PTX3.

UniProt ID P26022

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)