

PTX3 monoclonal antibody (MNB4)

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 18

[View Online »](#)

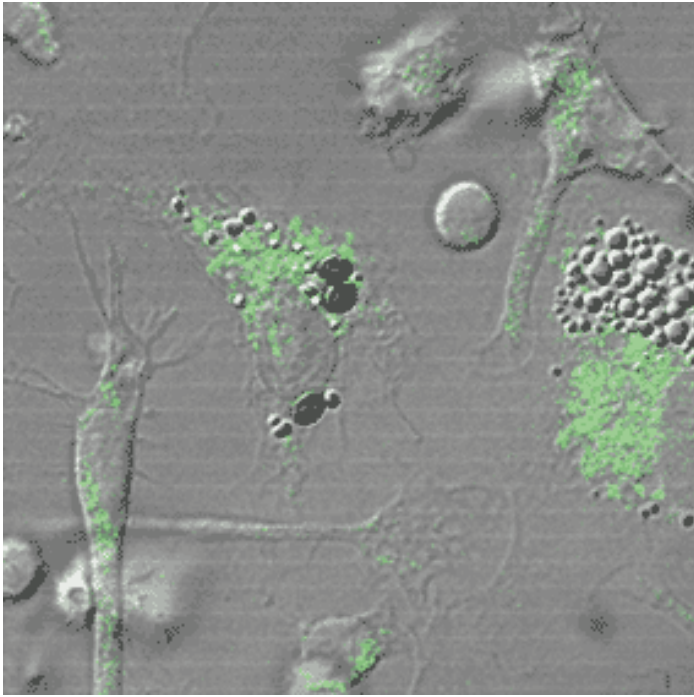
Ordering Information

[Order Online »](#)

ALX-804-464-C100	100µg
------------------	-------

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, TSG-14, TNF stimulated gene-14
Application	ELISA, Flow Cytometry, IHC (FS), IP, WB
Application Notes	Excellent for ELISA (capture).
Clone	MNB4
Formulation	Liquid. In PBS containing 0.1% sodium azide.
Host	Rat
Immunogen	Recombinant human PTX3 (pentraxin 3).
Isotype	IgG2a
Recommendation Dilutions/Conditions	ELISA (capture)Immunohistochemistry (frozen sections, 1µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Rat
Specificity	Recognizes the N-terminus of PTX3.
UniProt ID	P26022

Worry-free Guarantee This antibody is covered by our [Worry-Free Guarantee](#).



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)