PTX3 monoclonal antibody (MNB1)

This antibody is covered by our Worry-Free Guarantee.

Citations: 6

View Online »

Ordering Information

Order Online »

ALX-804-463-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 IHC (FS), IHC (PS), WB

Application Notes Excellent for Western blot.

Clone MNB1

Formulation Liquid. Purified antibody in PBS containing 0.1% sodium

azide.

Host Rat

Immunogen Recombinant human PTX3 (pentraxin 3).

lsotype lgG2b

Recommendation Dilutions/Conditions Western Blot (1µg/ml)Suggested dilutions/conditions may

not be available for all applications. Optimal conditions must be determined individually for each application.

Species Reactivity Human, Mouse

Specificity Recognizes the C-terminus of PTX3.

UniProt ID P26022

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

•

Last modified: May 29, 2024



info-

eu@enzolifesciences.com