

PTX3 polyclonal antibody (biotin conjugate)

Pentraxin 3 (PTX3), a member of the pentraxin family, is referred to as a long pentraxin and contains a unique PTX3 domain not found in C-reactive protein (CRP) or (Serum Amyloid P component) SAP. In contrast to CRP, PTX3 is produced from the major cell types involved in atherosclerotic lesions, namely vascular endothelial cells, vascular smooth muscle cells, macrophages, and neutrophils in response to inflammatory stimuli. Moreover, CRP is produced by the liver and represents a systemic response to local inflammation, whereas PTX3 is rapidly produced directly from damaged tissues and directly reflects the inflammatory state of the vasculature. PTX3 levels have been reported to be significantly elevated in acute myocardial infarction, and accordingly, statin treatment significantly decreases PTX3 levels. Due to the fact that PTX3 is able to reflect acute coronary syndrome (ACS) condition better than CRP, it is highly possible that PTX3 is a superior biomarker to predict future cardiovascular events.

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

Citations: 10

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-210-365B-C050	Biotin - 50µg
-------------------	---------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Once thawed store at +4°C and do not re-freeze.
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, ICC, IHC
Application Notes	Excellent for use as detection antibody in an ELISA format together with mAb to PTX3 (Prod. No. ALX-804-464 (capture)) and PTX3 standard (Prod. No. ALX-201-124).
Crossreactivity	Weakly cross-reacts with bovine PTX3.
Formulation	Liquid. In PBS containing 0.1% sodium azide.
Host	Rabbit
Immunogen	Recombinant human PTX3.
Purity Detail	Affinity purified.
Recommendation Dilutions/Conditions	ELISA (detection, 1:20,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Chicken, Human, Mouse, Rat
Specificity	Recognizes human, mouse and rat PTX3 and chicken homolog PPTX.
UniProt ID	P26022



ENZO LIFE SCIENCES,
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 1 46 60 42 00
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com

This antibody is covered by our Worry Free Guarantee