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 nutrophils 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

Last modified: May 29, 2024



info-

eu@enzolifesciences.com