

PTX3 (human) detection set [for ELISA application]

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.

Citations: 21

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Once thawed store biotinylated antibody 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, TSG14, TNF stimulated Gene-14
Application	ELISA
Application Notes	For the quantitative determination of PTX3 in human and bovine (weak) cell culture media, EDTA-plasma or serum (weak) samples.
Contents	Contains enough material to prepare 10 plates: – 1 vial (10µg) PTX3 (human) (rec.) (Prod. No. ALX-201-124) – 1 vial (50µg) PTX3, pAb (Biotin) [Detection Antibody] (Prod. No. ALX-210-365B) – 1 vial (100µg) PTX3, mAb (MNB4) [Capture Antibody] (Prod. No. ALX-804-464)
Sensitivity	Detection limit: 75pg/ml.
UniProt ID	P26022

Last modified: May 29, 2024



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

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

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

France
Phone: +33 472 440 655
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