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 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.

Citations: 21

View Online »

Ordering Information

Order Online »

ALX-850-299-KI01

1Set

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

