

TXB2 Polyclonal Antibody

(Catalog No. A-0400)

Background:

Thromboxane B2 (TXB2) is an inactive metabolite/product of thromboxane A2 (TXA2) - an enzyme involved in platelet aggregation, vasoconstriction and reproductive functions, but also known to be highly unstable under physiological conditions due to its very short half-life. TXB2 and its metabolite 2,3-dinor-TXB2 are released in abundant quantities during platelet aggregation and are commonly used to indirectly measure TXA2 production *in vivo*. A major metabolite of TXB2, 11-dehydro-TXB2, found in urine and plasma is produced via dehydrogenation by the enzyme 11-OH-dehydrogenase.

Description

Sheep polyclonal antibody

Formulation

Liquid. Contains sodium azide.

Immunogen

TXB2 conjugated to bovine serum albumin.

Specificity

Species Independent

Purification

Serum

Storage

Store at -20°C. Avoid multiple freeze/thaw cycles.

Alternative Names

Thromboxane B2

Application

ELISA

Ordering Information

Product	Size	Cat. No.
TXB2 Polyclonal Antibody	100 µl	A-0400-100

This product is for research purposes only. Not intended for use in diagnostic procedures.