## TXB<sub>2</sub> ELISA kit

Highly sensitive, well-validated ELISA kit with results in <3h

The TXB $_2$  ELISA kit is a colorimetric competitive enzyme immunoassay kit with results in < 3 hours. Absorbance is read at 405 nm. Detect COX I inhibition by measuring the production of downstream TXB $_2$  in a convenient 96 well plate based assay. This kit is highly cited in peer-reviewed publications in a diverse assortment of sample types including mucosa, tissue, serum, plasma, urine, and much more. The TXB $_2$  ELISA kit provides easy-to-use stable liquid color-coded reagents that reduce errors.

Thromboxane  $A_2$  (TXA $_2$ ) is involved in platelet aggregation, vasoconstriction and reproductive functions, but has a half life of only 37 seconds under physiological conditions. Thromboxane  $B_2$  (TXB $_2$ ) is the stable product of the non-enzymatic hydration of TXA2, thus the production of TXA $_2$  *in vivo* is typically monitored by measurement of TXB $_2$  and its metabolite 2,3-dinor TXB2. 11-dehydro-TXB2 is a major metabolite of Thromboxane  $B_2$  (TXB $_2$ ) found in urine and plasma produced by its dehydrogenation by the enzyme 11-OH-dehydrogenase.

Citations: 55

View Online »

**Ordering Information** 

Order Online »

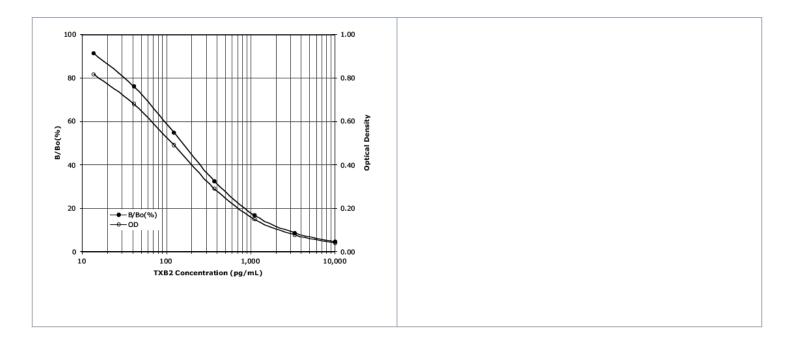
ADI-900-002

96 wells

Manuals, SDS & CofA

View Online »

- High sensitivity measurement, detecting as little as 10.54 pg/ml of TXB<sub>2</sub>
- High throughput format with results in < 3 hours for up to 37 samples in duplicate
- Widely cited in peer-reviewed publications in a diverse assortment of sample types including mucosa, tissue, serum, plasma, urine, and much more
- Reliable and reproducible results day-after-day and lot-after-lot



## **Handling & Storage**

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

**Product Details** 

Alternative Name Thromboxane B2

**Application** Colorimetric detection, ELISA

**Application Notes** For the quantitative determination of TXB<sub>2</sub> in culture

supernatants, plasma, serum, saliva, and urine from any species. Cited sample types include cell lysate, coronary effluent, liver perfusate, platelets, and whole blood.

Assay Time <3 hours

Compatibility This product is compatible with the Absorbance 96 Plate

Reader.

**Contents** GxR IgG Microtiter plate, Conjugate, Antibody, Assay

buffer, Wash buffer concentrate, Standard, pNpp

Substrate, Stop solution

**Sensitivity** 10.54 pg/ml (range 13.7 – 10,000 pg/ml)

Species Reactivity Species independent

Wavelength 405 nm

Last modified: May 29, 2024

