## TXB2 ELISA kit

## The TXB<sub>2</sub> ELISA kit is a colorimetric competitive enzyme immunoassay kit

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** »

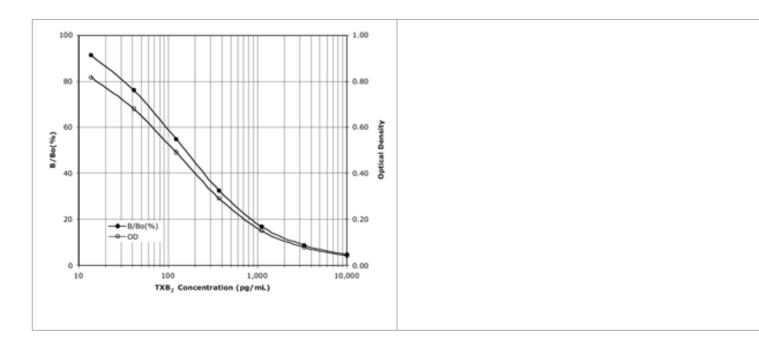
ADI-901-002

5x96 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
- Save with bulk kit package!



## **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.

Compatibility This product is compatible with the Absorbance 96 Plate

Reader.

Contents GxR Microtiter plate, Conjugate, Antibody, Assay buffer

concentrate, 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

