

TXB₂ ELISA kit

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

The TXB₂ 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 TXB₂ 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 TXB₂ ELISA kit provides easy-to-use stable liquid color-coded reagents that reduce errors.

Thromboxane A₂ (TXA₂) is involved in platelet aggregation, vasoconstriction and reproductive functions, but has a half life of only 37 seconds under physiological conditions. Thromboxane B₂ (TXB₂) is the stable product of the non-enzymatic hydration of TXA₂, thus the production of TXA₂ *in vivo* is typically monitored by measurement of TXB₂ and its metabolite 2,3-dinor TXB₂. 11-dehydro-TXB₂ is a major metabolite of Thromboxane B₂ (TXB₂) found in urine and plasma produced by its dehydrogenation by the enzyme 11-OH-dehydrogenase.

Citations: 56

[View Online »](#)

Ordering Information

[Order Online »](#)

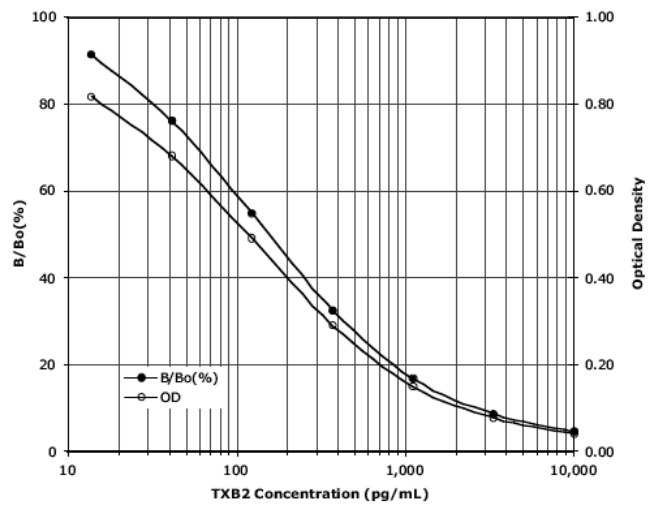
ADI-900-002

96 wells

Manuals, SDS & CofA

[View Online »](#)

- High sensitivity measurement, detecting as little as 10.54 pg/ml of TXB₂
- 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 TXB₂ 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 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



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