

TXB₂ ELISA kit

The TXB₂ ELISA kit is a colorimetric competitive enzyme immunoassay kit

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: 51

[View Online »](#)

Ordering Information

[Order Online »](#)

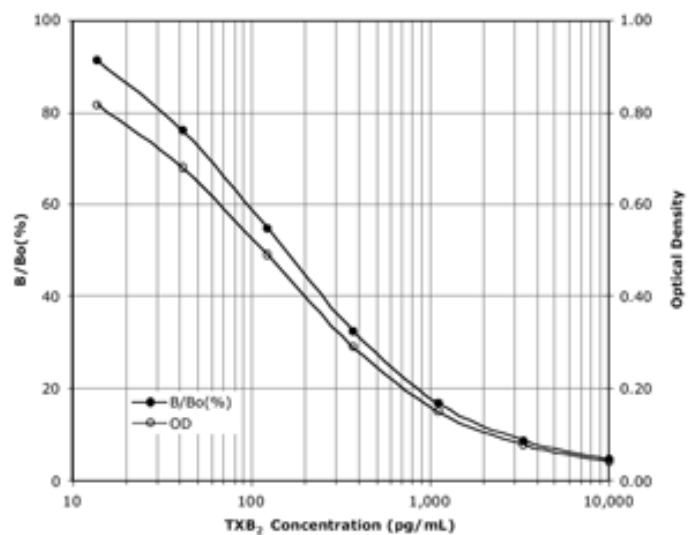
ADI-901-002

5x96 wells

Manuals, SDS & CofA

[View Online »](#)

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



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)