

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: 52

[View Online »](#)

Ordering Information

[Order Online »](#)

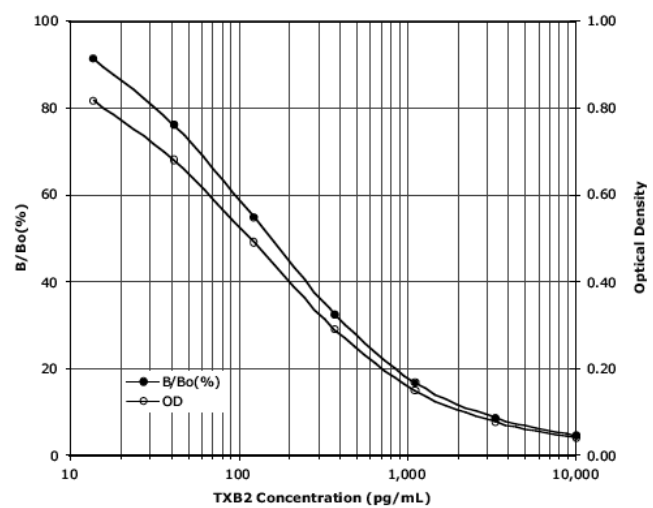
ADI-900-002

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.

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



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)