

11-dehydro-TXB2

ELISA kit

The 11-dehydro-TXB₂ ELISA kit is a colorimetric competitive enzyme immunoassay kit

The 11-dehydro-TXB₂ EIA kit is a colorimetric competitive enzyme immunoassay kit with results in 3 hours. Absorbance is read at 405 nm. This kit offers ultrasensitive quantification of 11-d-TXB₂ in a variety of sample matrices with negligible reactivity with related molecules in a convenient 96-well format offers higher throughput analysis compared to conventional analytical methods.

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: 10

[View Online »](#)

Ordering Information

[Order Online »](#)

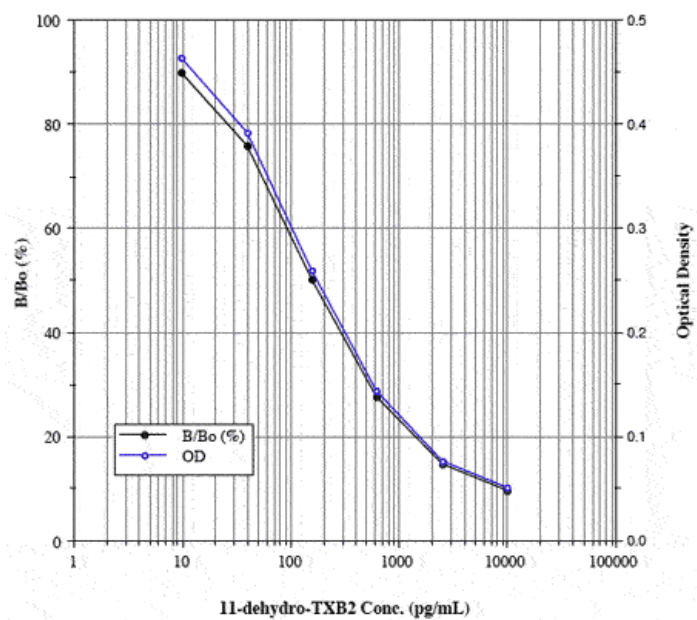
ADI-900-092

96 wells

Manuals, SDS & CofA

[View Online »](#)

- Higher throughput - measure up to 38 samples in duplicate in just 3 hours
- Easy-to-use - works with any standard colorimetric plate reader, no specialized training required
- Ready-to-use liquid color-coded reagents save time and reduce errors
- More sensitive than other commercially available kits
- Specific - low reactivity to structurally related molecules



Handling & Storage

Use/Stability Store all components at +4°C, except standard and conjugate at -20°C.

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 11-dehydro-Thromboxane B2

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of 11-dehydro-TXB₂ in culture supernatants and urine from any species. Cited samples include serum.

Assay Time 3 hours

Compatibility This product is compatible with the [Absorbance 96 Plate Reader](#).

Contents GxR IgG Microtiter plate, Assay buffer, Conjugate, Antibody, Wash buffer concentrate, Standard, pNpp Substrate, Stop solution

Sensitivity 4.31 pg/ml (range 9.8 – 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-fr@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