

# 11-dehydro-TXB2

## ELISA kit

The 11-dehydro-TXB<sub>2</sub> ELISA kit is a colorimetric competitive enzyme immunoassay kit

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

Citations: 10

[View Online »](#)

### Ordering Information

[Order Online »](#)

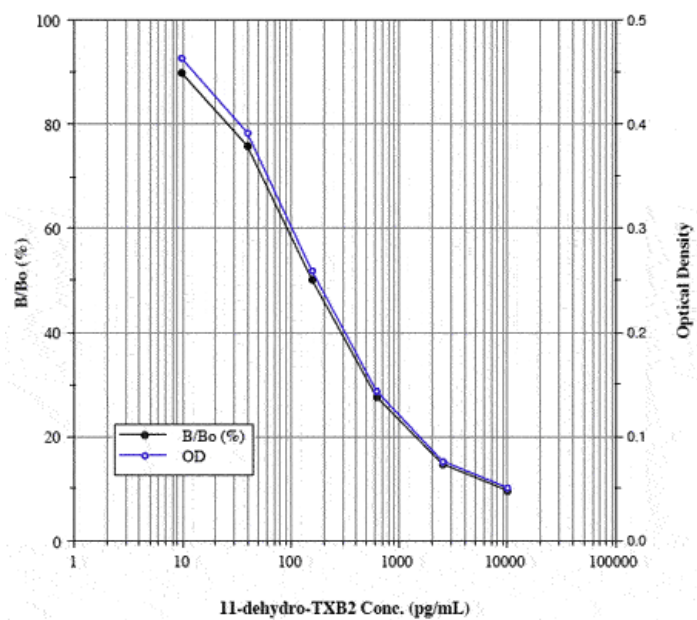
ADI-900-092

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-TXB<sub>2</sub> 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



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-  
be@enzolifesciences.com](mailto:info-be@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)