

Product datasheet

8 isoprostane ELISA Kit ab175819

[1 References](#) [1 Image](#)

Overview

Product name 8 isoprostane ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Overall	12			0.43%

Inter-assay

Sample	n	Mean	SD	CV%
Overall	12			1.75%

Sample type Urine, Serum, Plasma, Cell culture extracts, Tissue

Assay type Competitive

Sensitivity 1 pg/ml

Range 0.005 ng/ml - 5 ng/ml

Assay duration Multiple steps standard assay

Product overview

Abcam's 8 isoprostane competitive *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the determination of 8 isoprostane levels in biological samples. The isoprostanes are a family of eicosanoids of non-enzymatic origin produced by the random oxidation of tissues phospholipids by oxygen radicals. A recent NIH-sponsored study on Biomarkers of Oxidative Stress has indicated that 8 isoprostane is the best index of oxidative injury in a well-accepted oxidant stress rat model. In addition, plasma 8 isoprostane levels were found to be elevated in elderly subjects with severe hypertension and in urine from subjects with high fat diet-induced liver steatosis.

This kit can be used for the determination of 8 isoprostane in diluted urine, serum, plasma, cells and tissues following proper isolation and purification of the eicosanoid from the isoprostane-containing sample.

Tested applications **Suitable for:** Competitive ELISA

Platform Microplate

Properties

Storage instructions

Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
8-isoprostane HRP Conjugates	1 x 12µl
8-isoprostane Standard	1 x 2µl
HRP Buffer	1 x 15ml
Pre-coated 96-Well Microplate	1 unit
Sample Dilution Buffer	1 x 25ml
TMB Substrate	1 x 24ml
Wash Buffer Solution	1 x 25ml

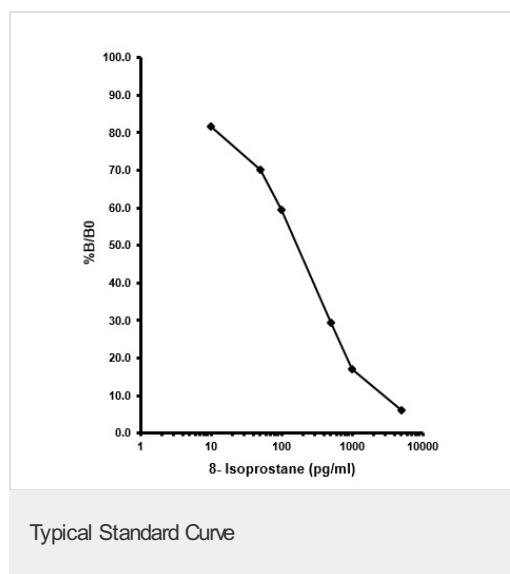
Applications

Our [Abpromise guarantee](#) covers the use of **ab175819** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
Competitive ELISA		Use at an assay dependent concentration.

Images



Example of standard curve obtained with ab175819.

Typical Standard Curve

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours

- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <http://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors