

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

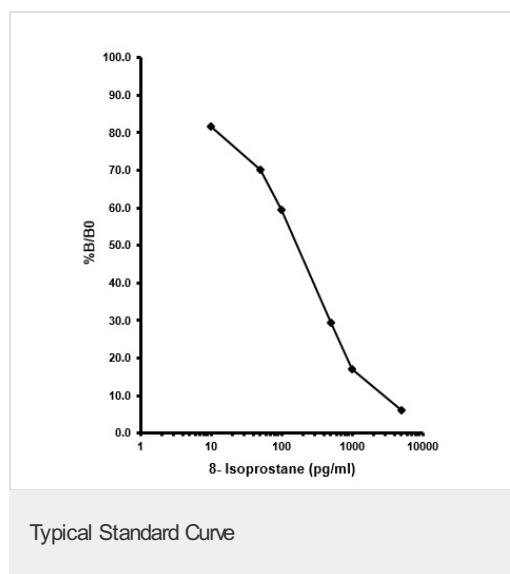
Applications

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**.

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 <https://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