

## Product datasheet

# 14,15 DHET ELISA Kit ab175811

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

### Overview

**Product name** 14,15 DHET ELISA Kit

**Detection method** Colorimetric

**Precision**

Intra-assay

Sample	n	Mean	SD	CV%
Overall	12			0.93%

Inter-assay

Sample	n	Mean	SD	CV%
Overall	12			1.31%

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

**Assay type** Competitive

**Sensitivity** = 3 pg/ml

**Range** 0.1 ng/ml - 1000 ng/ml

**Assay duration** Multiple steps standard assay

**Product overview** Abcam's 14,15 DHET competitive *in vitro* ELISA Kit is designed for accurate quantitative measurement in serum, plasma, urine, cell culture, tissue samples following proper isolation and purification.

This competitive ELISA kit is based on competition between the 14,15-DHET epitope and the 14,15-DHET-HRP conjugate for a limited number of binding sites available from the anti-14,15-DHET antibody; which is coated to the wells of the 96 well ELISA plate. The conjugate concentration is held as a constant in each well, while the concentration of the 14,15-DHET is variable, based on the concentration of the sample or standard. Thus the amount of the 14,15-DHET conjugate which is able to bind to each of the wells is inversely proportional to the concentration of 14,15-DHET in the standard or sample. The amount of the conjugate which is bound to each well is then determined by the amount of color obtained, when TMB is added. The TMB reacts with the HRP available in the well. With the addition of sulfuric acid, the blue colored product is converted into a yellow colored product, which can be read on a plate reader at 450 nm.

This competitive ELISA kit is for determination of 14,15-DHET levels in biological samples. Specificity of the ELISA has been published. The 14,15-DHET level exhibited strong positive correlation with hypertension, brain injury and stroke in rodents. 14,15-DHET is a representative metabolite of soluble epoxide hydrolase-mediated metabolism of EETs, which are generated by arachidonic acid epoxygenase activity of cytochromes P450. Human blood 14,15-DHET levels were measured using the 14,15-DHET ELISA kit. Increased 14,15-DHET levels of human cells as detected by the 14,15-DHET ELISA were indicative of the neoplastic and metastatic phenotype of carcinoma cells.

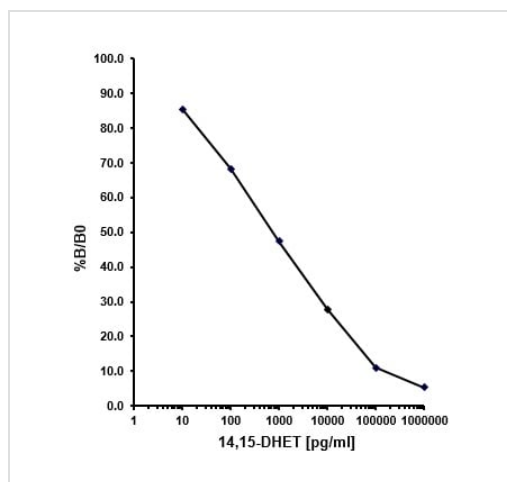
**Platform** Microplate

## Properties

**Storage instructions** Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
14,15-DHET Standard	1 x 2µl
14,15-DHET-HRP Conjugates	1 x 12µl
HRP Buffer	1 x 15ml
Pre-coated 96-Well Microplate	1 unit
10X Sample Dilution Buffer	1 x 25ml
TMB Substrate	1 x 24ml
10X Wash Buffer Solution	1 x 25ml

## Images



Representative standard curve using ab175811 14,15 DHET ELISA Kit.

Typical Standard Curve

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

### **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