

## Product datasheet

# Human FABP2 ELISA Kit ab193700

★★★★★ 1 Abreviews 1 Image

### Overview

**Product name** Human FABP2 ELISA Kit

**Detection method** Colorimetric

**Precision**

Intra-assay

Sample	n	Mean	SD	CV%
--------	---	------	----	-----

Overall < 10%

Inter-assay

Sample	n	Mean	SD	CV%
--------	---	------	----	-----

Overall < 12%

**Sample type** Serum, Plasma, Cell culture media

**Assay type** Sandwich

**Sensitivity** 25 pg/ml

**Range** 25 pg/ml - 100000 pg/ml

**Recovery**

Sample specific recovery

Sample type	Average %	Range
-------------	-----------	-------

Serum 129.8 120% - 136%

Plasma 126.5 110% - 135%

Cell culture media 112.4 103% - 123%

**Assay duration** Multiple steps standard assay

**Species reactivity** **Reacts with:** Human

**Product overview**

Abcam's FABP2 Human ELISA Kit (ab193700) is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of human FABP2 in sera, plasma and cell culture supernatants.

This assay employs an antibody specific for human FABP2 coated on a 96-well plate. Standards

and samples are pipetted into the wells and the immobilized antibody captures FABP2 present in the samples. The wells are washed and biotinylated anti-human FABP2 antibody is added. After washing away any unbound biotinylated antibody, an HRP-conjugated streptavidin is pipetted to the wells. After incubation, the wells are again washed, followed by the addition of a TMB substrate solution to the wells. Color will develop in proportion to the amount of FABP2 bound in each well. Addition of the Stop Solution will change the color from blue to yellow, and the intensity of the color is measured at 450 nm.

**Platform** Pre-coated microplate (12 x 8 well strips)

## Properties

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

Components	1 x 96 tests
1500X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer Concentrate	1 x 25ml
5X Assay Diluent B	1 x 15ml
Assay Diluent C	1 x 30ml
Biotinylated Human FABP2 detection antibody (lyophilized)	2 vials
Human FABP2 Standards (lyophilized)	2 vials
Pre-coated Human FABP2 Microplate (12 strips x 8 wells)	1 unit
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

**Function** FABP are thought to play a role in the intracellular transport of long-chain fatty acids and their acyl-CoA esters. FABP2 is probably involved in triglyceride-rich lipoprotein synthesis. Binds saturated long-chain fatty acids with a high affinity, but binds with a lower affinity to unsaturated long-chain fatty acids. FABP2 may also help maintain energy homeostasis by functioning as a lipid sensor.

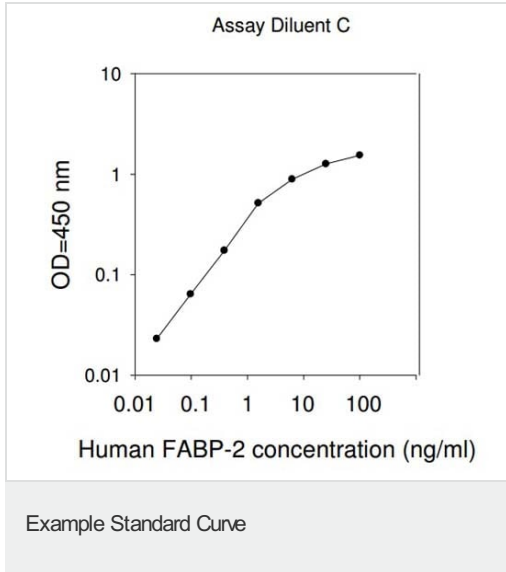
**Tissue specificity** Expressed in the small intestine and at much lower levels in the large intestine. Highest expression levels in the jejunum.

**Sequence similarities** Belongs to the calycin superfamily. Fatty-acid binding protein (FABP) family.

**Domain** Forms a beta-barrel structure that accommodates the hydrophobic ligand in its interior.

**Cellular localization** Cytoplasm.

## Images



**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