

### Rat L-FABP ELISA Kit ab190545

1 Image

#### Overview

**Product name** Rat L-FABP ELISA Kit

**Detection method** Colorimetric

**Precision**

Intra-assay

Sample	n	Mean	SD	CV%
Overall				< 10%

Inter-assay

Sample	n	Mean	SD	CV%
Overall				< 10%

**Sample type** Serum, Plasma

**Assay type** Sandwich (quantitative)

**Sensitivity** 1.278 ng/ml

**Range** 3.125 ng/ml - 200 ng/ml

**Recovery**

Sample specific recovery

Sample type	Average %	Range
Serum	> 85	% - %

**Assay duration** Multiple steps standard assay

**Species reactivity** **Reacts with:** Rat

**Product overview**

Abcam's L-FABP Rat ELISA Kit (ab190545) is an *in vitro* enzyme-linked immunosorbent assay (ELISA) for the quantitative measurement of L-FABP in rat serum and plasma.

In this assay the L-FABP present in samples reacts with the anti-L-FABP antibodies which have been adsorbed to the surface of polystyrene microtitre wells. After the removal of unbound proteins by washing, anti-L-FABP antibodies conjugated with horseradish peroxidase (HRP) are added. These enzyme-labeled antibodies form complexes with the previously bound L-FABP. Following another washing step, the enzyme bound to the immunosorbent is assayed by the addition of a chromogenic substrate, 3,3',5,5'-tetramethylbenzidine (TMB). The quantity of bound enzyme varies directly with the concentration of L-FABP in the sample tested; thus, the

absorbance, at 450 nm, is a measure of the concentration of L-FABP in the test sample. The quantity of L-FABP in the test sample can be interpolated from the standard curve constructed from the standards, and corrected for sample dilution.

**Platform** Microplate

### Properties

Components	1 x 96 tests
1X Diluent Concentrate	1 x 60ml
20X Wash Buffer Concentrate	1 x 50ml
Chromogen Substrate Solution	1 x 12ml
L-FABP Rat Antibody coated microwells	1 x 96 tests
L-FABP Rat Calibrator (Lyophilized)	1 vial
L-FABP Rat HRP Conjugate	1 vial
Stop Solution	1 x 12ml

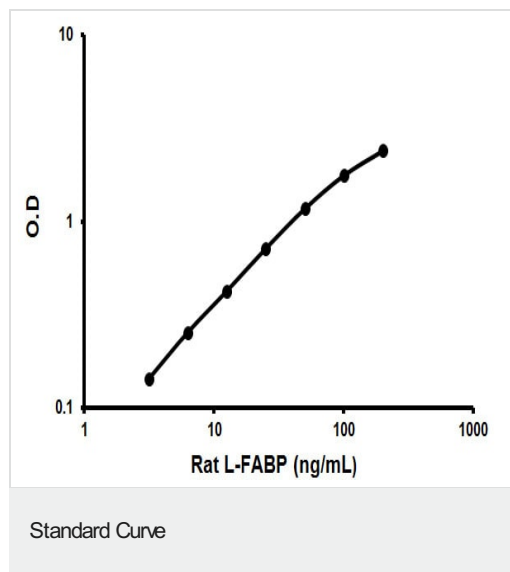
**Function** Binds free fatty acids and their coenzyme A derivatives, bilirubin, and some other small molecules in the cytoplasm. May be involved in intracellular lipid transport.

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

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

**Cellular localization** Cytoplasm.

### Images



Representative standard curve using ab190545 L-FABP Rat ELISA Kit.

**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