

Human Fatty Acid Oxidation In-Cell ELISA Kit **ab118182**

[3 References](#) [4 Images](#)

Overview

Product name	Human Fatty Acid Oxidation In-Cell ELISA Kit
Detection method	Colorimetric
Sample type	Adherent cells, Suspension cells
Assay type	Cell-based (quantitative)
Assay duration	Multiple steps standard assay
Product overview	ab118182 uses quantitative immunocytochemistry to measure protein levels or post-translational modifications in cultured cells. Cells are fixed in a microplate and targets of interest are detected with highly specific, well-characterized monoclonal antibodies, and levels are quantified with IRDye®-labeled Secondary Antibodies. IR imaging and quantitation is performed using a LI-COR® Odyssey® or Aeries® system.

Plates are available in our ICE (In-Cell ELISA) Support Pack (**ab111542**) which can be bought separately.

Notes Upon receipt spin down the contents of the IRDye®-labeled Secondary Antibody tube and protect from light. Store all components upright at 4°C. This kit is stable for at least 6 months from receipt.

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances.

It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

Platform Microplate

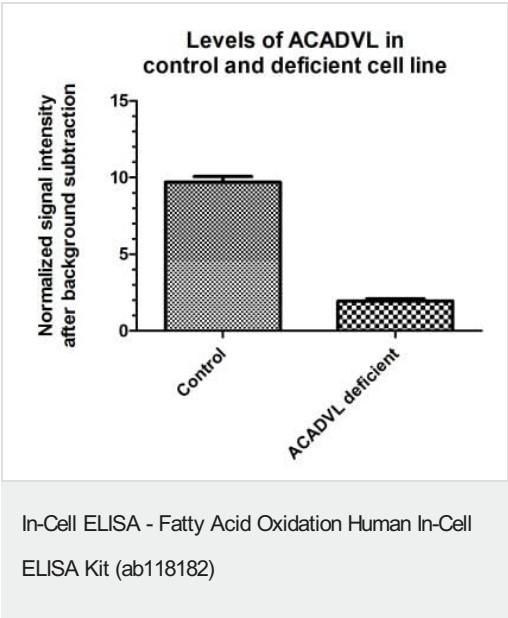
Properties

Storage instructions Store at +4°C. Please refer to protocols.

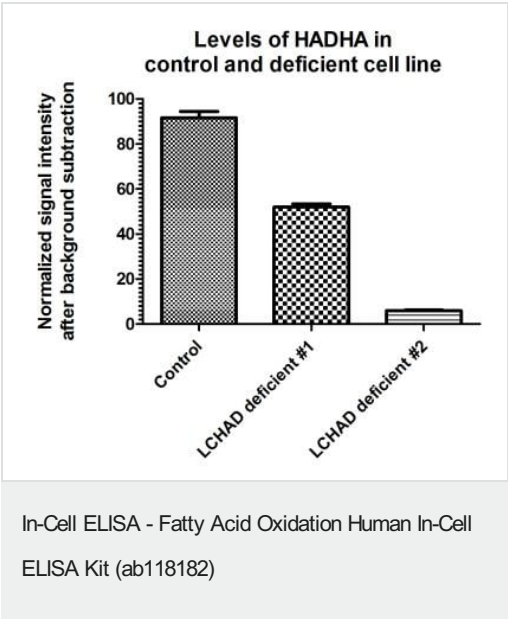
Components	1 x 96 tests
1000X IRDye 800-labeled Secondary antibody	1 x 0.04ml
100X ACADM Primary Antibody	1 x 0.12ml
100X ACADVL Primary Antibody	1 x 0.12ml

Components	1 x 96 tests
100X HADHA Primary Antibody	1 x 0.12ml
10X Blocking Buffer	1 x 20ml
10X Phosphate Buffered Saline	1 x 100ml
33x Triton X-100	1 x 5ml
400X Tween-20	1 x 4ml
1X Janus Green Stain	1 x 17ml

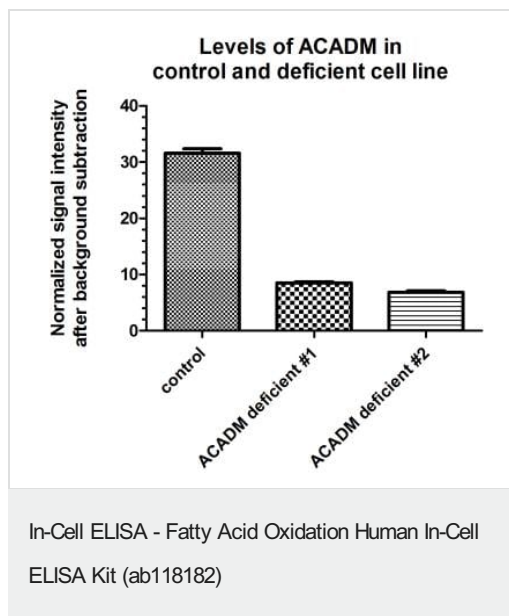
Images



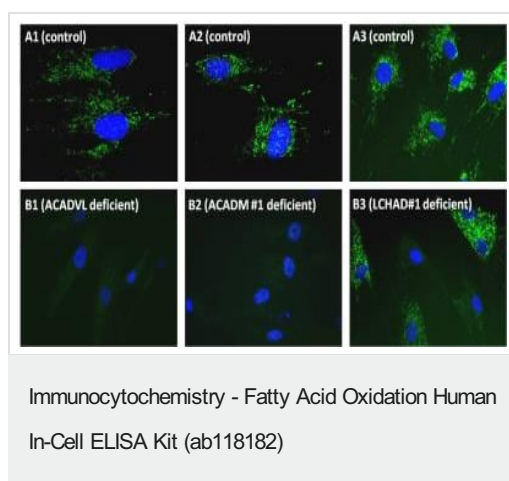
Panel A shows levels of ACADVL protein in a deficient cell line (ACADVL:p.[N122D]).



Panel C shows levels of HADHA protein in two deficient cell lines: (1) HADHA:p.[E474Q] and (2) HADHB:p. [R61H];[R247H]. Note that although LCHAD#1 deficient has a partial deficiency as observed by ICC.



Panel B shows levels of ACADM protein in two deficient cell lines (ACADM:p. [K304E]).



Antibody specificity demonstrated by immunocytochemistry. Visualization under the microscope was carried out with a completely opened aperture and a very narrow field of visualization at 40x. Panel A shows control fibroblasts and panel B shows deficient fibroblasts. Left panel shows staining with anti-ACADVL ab, center panel with anti-ACADM ab and right panel with anti-HADHA ab. Note that although LCHAD#1 deficient had a characterized homozygous mutation, the antibody shows a mosaic pattern of staining with about 1/3 of the cells lacking HADHA staining and an overall 40% reduction of signal as observed by ICE.

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