

Fatty Acid Oxidation Assay Kit (flow cytometry) ab118183

9 References 7 图像

概述

产品名称	Fatty Acid Oxidation Assay试剂盒(flow cytometry)
检测方法	Fluorescent
样品类型	Adherent cells, Suspension cells
检测类型	Quantitative
产品概述	Fatty Acid Oxidation Assay Kit ab118183 contains antibodies against key enzymes of the mitochondrial fatty acid oxidation pathway for flow cytometry. The assay combines the power of single cell analysis obtained from flow cytometry with the specificity of antibody-based immunostaining to quantify protein levels in cultured cells. Cells are harvested and fixed/permeabilized in suspension, targets of interest are detected indirectly with highly specific well-characterized monoclonal antibodies that are then labeled with fluorescent antibodies

说明	<p>This product was previously called Fatty Acid Oxidation Human Flow Cytometry Kit.</p> <p>Spin down the contents of the antibody vials upon receipt of the kit. Store all components upright at 4°C. This kit is stable for at least 6 months from receipt.</p> <p>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.</p> <p>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.</p>
----	--

平台	Flow cytometer
----	----------------

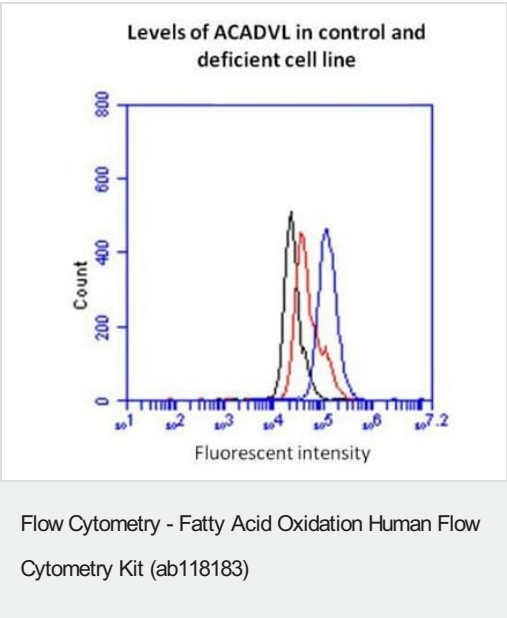
性能

存放说明	Store at +4°C. Please refer to protocols.
------	---

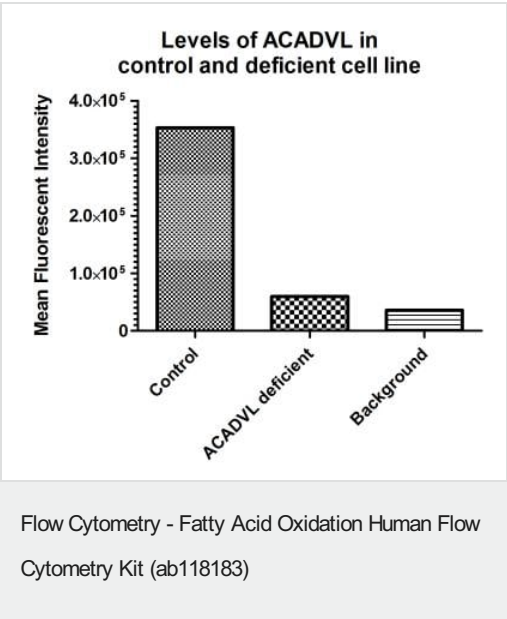
组件	96 tests
100X ACADM Primary Antibody	1 x 0.12ml
100X ACADVL Primary Antibody	1 x 0.12ml

组件	96 tests
100X HADHA Primary Antibody	1 x 0.12ml
100X Normal Mouse IgG	1 x 0.12ml
100X Triton X-100	1 x 1.25ml
10X Phosphate Buffered Saline	1 x 100ml
400X Tween-20	1 x 4ml
Blocking Solution	1 x 25ml

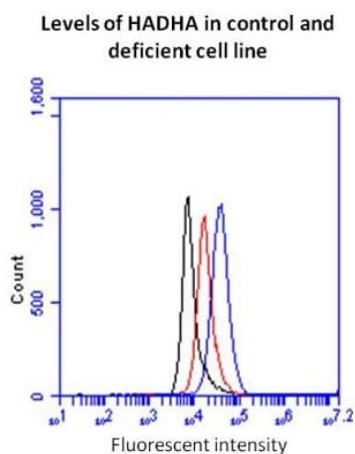
图片



Expression levels of ACADVL was determined in primary fibroblasts by flow cytometry. Blue: control cell line. Red: deficient cell lines extracted from patients with well characterized mutations in each of the enzymes ACADVL:p[N122D]. Black: Background fluorescence was determined with the supplied negative control.

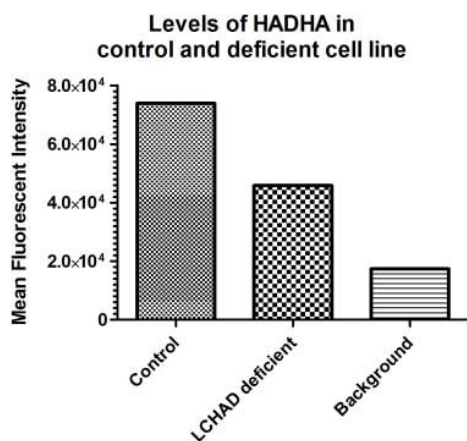


After background subtraction, the ACADVL deficient cell line shows a 93% reduction in the level of the ACADVL protein.



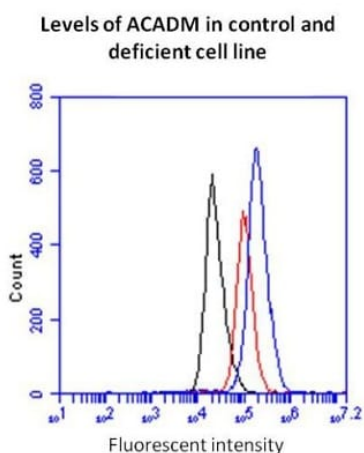
Flow Cytometry - Fatty Acid Oxidation Human Flow Cytometry Kit (ab118183)

Expression levels of HADHA was determined in primary fibroblasts by flow cytometry. Blue: control cell line. Red: deficient cell lines extracted from patients with well characterized mutations in each of the enzymes HADHB:p[R61H];[R247H]. Black: Background fluorescence was determined with the supplied negative control.



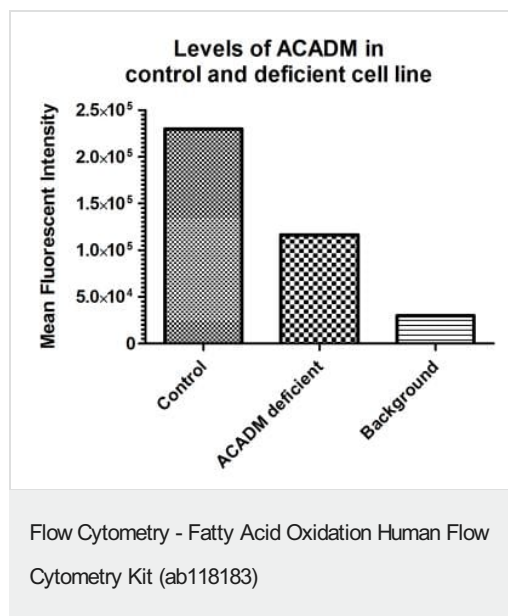
Flow Cytometry - Fatty Acid Oxidation Human Flow Cytometry Kit (ab118183)

After background subtraction, the LCHAD deficient cell line shows a 50% reduction in the level of the HADHA protein.

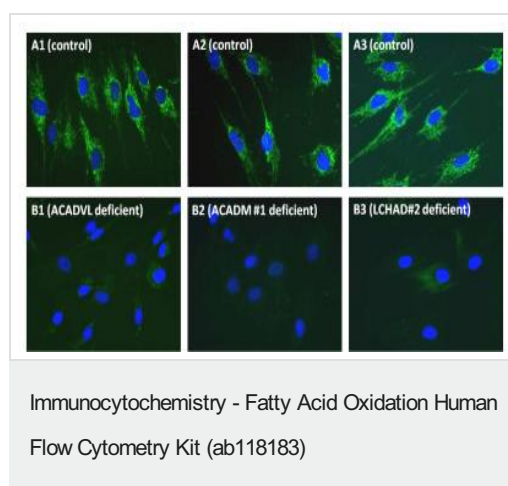


Flow Cytometry - Fatty Acid Oxidation Human Flow Cytometry Kit (ab118183)

Expression levels of ACADM was determined in primary fibroblasts by flow cytometry. Blue: control cell line. Red: deficient cell lines extracted from patients with well characterized mutations in each of the enzymes ACADM:p[K604E] Black: Background fluorescence was determined with the supplied negative control.



After background subtraction, the ACADM deficient cell line shows a 57% reduction in the level of the ACADM protein.



Antibody specificity demonstrated by immunocytochemistry. Cells were processed with the Flow cytometry protocol as explained above. 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.

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.cn/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors