

Lysyl Oxidase Activity Assay Kit (Fluorometric) ab112139

34 References 4 图像

概述

产品名称	Lysyl Oxidase Activity Assay试剂盒(Fluorometric)
检测方法	Fluorescent
样品类型	Cell culture supernatant, Cell culture extracts, Tissue Extracts
检测类型	Semi-quantitative
灵敏度	40 ng/well
检测时间	0h 30m
产品概述	<p>Lysyl Oxidase Activity Assay Kit / LOX Activity Assay Kit (Fluorometric) (ab112139) provides a simple method to measure lysyl oxidase (LOX) activity in cell and tissue extracts from mammals and other species, as well as physiological solutions.</p> <p>The LOX activity assay protocol uses a proprietary LOX substrate that releases hydrogen peroxide upon transformation by the LOX present in the sample. Hydrogen peroxide is in turn detected using a red fluorescence substrate for HRP-coupled reactions. This leads to increase in fluorescence that can be easily detected at Ex/Em = 540/590 nm in a fluorescence microplate reader.</p> <p>This assay is semi-quantitative as it does not contain a LOX standard for calibration. When a known concentration of LOX is used, the assay can detect activity from as low as 40 ng of lysyl oxidase in solution.</p> <p>The assay is highly sensitive and its unique detection method eliminates the interference that occurs in certain biological samples.</p>
说明	<p>Lysyl oxidase (protein-lysine-6-oxidase, LOX, EC 1.4.3.13) is an extracellular copper-dependent enzyme that catalyzes formation of aldehydes from lysine residues in collagen and elastin precursors.</p> <p>The activity of Lysyl oxidase in biological samples is traditionally assessed by tritium release end-point assays using radio isotope labeled collagen or elastin substrates.</p>
平台	Microplate reader

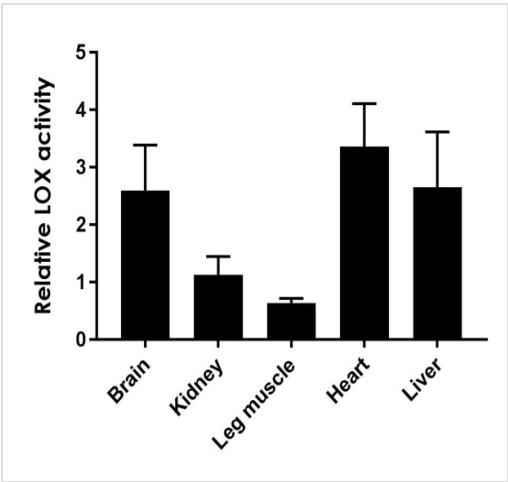
性能

存放说明

Store at -20°C. Please refer to protocols.

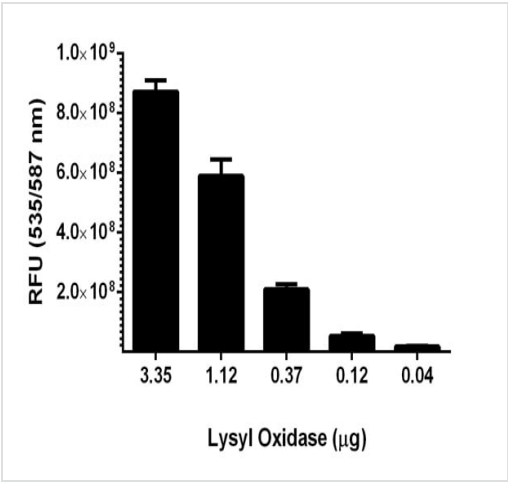
组件	500 tests
Assay Buffer	1 x 50ml
DMSO	1 x 200µl
Horseradish Peroxidase	1 vial
HRP Substrate	1 vial

图片



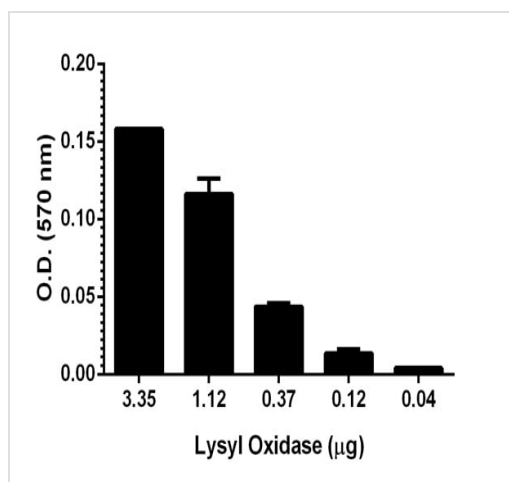
Relative LOX activity levels in mouse tissue. Tissue samples were prepared following assay protocol. Protein concentration was determined and samples were diluted 2-30 fold. LOX activity levels were measured after 15 minutes incubation in a fluorometric plate reader at Ex/Em = 535/587 nm. LOX activity levels are relative to background noise (blank control).

Lysyl Oxidase Activity Assay Kit (Fluorometric)
(ab112139)



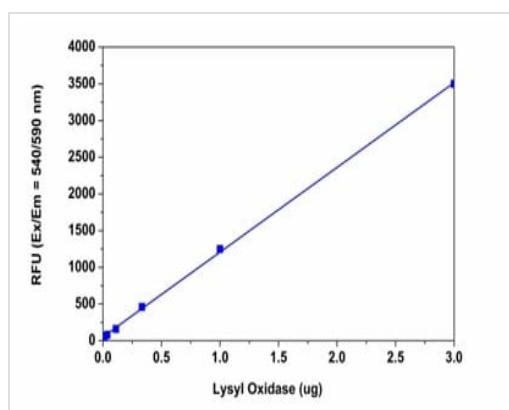
LOX activity fluorescence (RFU) values vs quantity (µg). Recombinant human LOX2 was serially diluted 3.35-0.04 (1/3) and activity was measured following assay procedure.

Functional Studies - Lysyl Oxidase Activity Assay
Kit (Fluorometric) (ab112139)



LOX activity absorbance (OD) values vs quantity (µg). Recombinant human LOX2 was serially diluted 3.35-0.04 (1/3) and activity was measured following assay procedure. Colorimetric detection is approximately 10-times lower than fluorometric.

Functional Studies - Lysyl Oxidase Activity Assay
Kit (Fluorometric) (ab112139)



Typical lysyl oxidase (LOX) dose response curve. Known amounts of LOX were added to wells and reaction was run following assay protocol. Fluorescence was measured on a solid black 96-well plate using a Gemini fluorescence microplate reader (Molecular Devices). Activity from as low as 40 ng/mL of LOX can be detected after 30 minutes incubation.

Lysyl Oxidase Activity Assay Kit (Fluorometric)
(ab112139)

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