


KDM1/LSD1 Activity Quantification Assay Kit (Colorimetric)

ab113459

2 References 2 图像

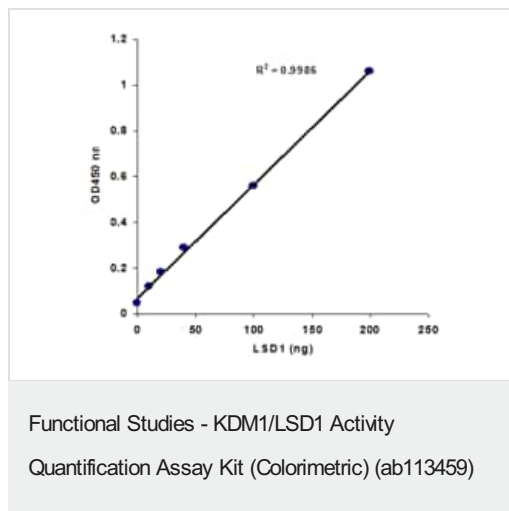
概述

产品名称	KDM1/LSD1 Activity Quantification Assay试剂盒(Colorimetric)
检测方法	Colorimetric
检测类型	Enzyme activity
检测时间	3h 00m
种属反应性	与反应: Mouse, Human, Plants 预测可用于: Mammals, Other species 
产品概述	<p>KMD1/LSD1 (Lysine Specific Histone Demethylase 1) is a histone demethylase that only acts on mono- or di-methylated lysines. KDM1/LSD1 acts as a co-repressor by mediating demethylation of H3(meK4), a specific tag for epigenetic transcriptional activation.</p> <p>KDM1/LSD1 Activity Quantification Assay Kit (Colorimetric) (ab113459) is designed for the easy and fast colorimetric measurement of KDM1/LSD1 activity or inhibition. The antibody-based, immunospecific method directly detects KDM1/LSD1-converted demethylated products, rather than by-products, in a 96 stripwell microplate format.</p> <p>This kit may cross react with LSD2.</p>
平台	Microplate reader

性能

存放说明 Please refer to protocols.

组件	48 tests	96 tests
10X Wash Buffer	1 x 14ml	1 x 28ml
8-Well Assay Strips (with Frame)	6 units	12 units
Adhesive Covering Film	1 unit	1 unit
Capture Antibody, 1000 µg/mL	1 x 5µl	1 x 10µl



Demonstration of high sensitivity of KDM1 / LSD1 activity assay achieved by using recombinant KDM1 / LSD1 with ab113459.

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