


KDM5/JARID Activity Quantification Assay Kit (Colorimetric) ab113463

2 References 1 Image

Overview

Product name	KDM5/JARID Activity Quantification Assay Kit (Colorimetric)
Detection method	Colorimetric
Assay type	Enzyme activity
Assay time	3h 00m
Species reactivity	Reacts with: Mouse, Human Predicted to work with: Mammals 
Product overview	<p>KDM5/JARID (Lysine-specific demethylase 5) is a family of histone demethyltransferases involved in H3(meK4) specific demethylation. KDM5/JARID demethylases are JmjC proteins that remove di- and tri-methylation from H3(meK4) but have no activity on mono-methylated residues.</p> <p>KDM5/JARID Activity Quantification Assay Kit (Colorimetric) (ab113463) is designed for the easy and fast colorimetric measurement of KDM5/JARID activity or inhibition. The antibody-based, immunospecific method directly detects KDM5/JARID-converted demethylated products, rather than by-products, in a 96 stripwell microplate format.</p>
Platform	Microplate reader

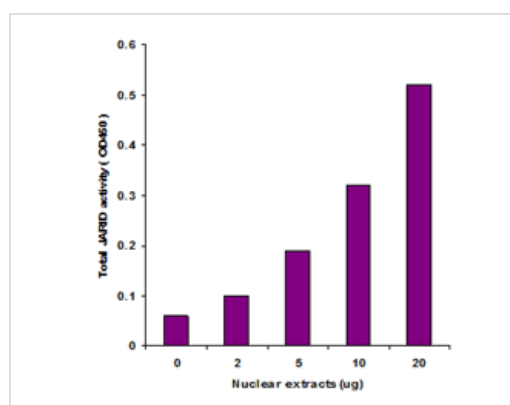
Properties

Storage instructions Please refer to protocols.

Components	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
Co-factor 1	1 x 30µl	1 x 60µl

Components	48 tests	96 tests
Co-factor 2	1 x 30µl	1 x 60µl
Co-factor 3	1 x 30µl	1 x 60µl
Detection Antibody, 400 µg/mL	1 x 6µl	1 x 12µl
Developer Solution	1 x 5ml	1 x 10ml
JARID Assay Buffer	1 x 4ml	1 x 8ml
JARID Assay Standard, 50 µg/mL	1 x 10µl	1 x 20µl
JARID Substrate, 50 µg/mL	1 x 60µl	1 x 120µl
Stop Solution	1 x 5ml	1 x 10ml

Images



Demonstration of high sensitivity of the KDM5/JARID activity assay achieved by using A549 nuclear extracts with ab113463.

Functional Studies - KDM5/JARID Activity
Quantification Assay Kit (Colorimetric) (ab113463)

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