

## Product datasheet

# KDM4/JMJD2 Activity Quantification Assay Kit (Fluorometric) ab113462

[1 References](#) [1 Image](#)

### Overview

<b>Product name</b>	KDM4/JMJD2 Activity Quantification Assay Kit (Fluorometric)
<b>Assay type</b>	Enzyme activity
<b>Assay time</b>	3h 00m
<b>Species reactivity</b>	<b>Reacts with:</b> Mouse, Human

**Predicted to work with:** Mammals

### Product overview

KDM4/JMJD2 (Lysine-specific demethylase 4) is a family of histone demethyltransferases involved in H3(meK9) specific demethylation. The KDM4/JMJD2 family are JmjC proteins that specifically removes tri-methylation from H3(meK9) and H3(meK36) but has no activity on mono- or di-methylated residues.

KDM4/JMJD2 Activity Quantification Assay Kit (Fluorometric) (ab113462) is designed for the easy and fast fluorometric measurement of KDM4/JMJD2 activity or inhibition. The antibody-based, immunospecific method directly detects KDM4/JMJD2-converted demethylated products, rather than by-products, in a 96 stripwell microplate format.

### Notes

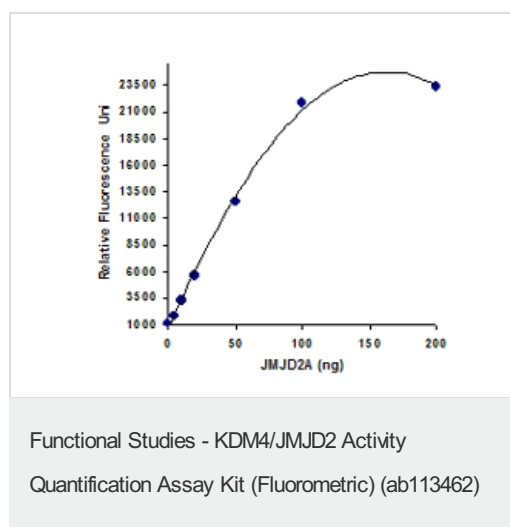
### Properties

**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
Fluoro Developer	1 x 10µl	1 x 20µl
Fluoro Dilution	1 x 4ml	1 x 8ml
Fluoro Enhancer	1 x 10µl	1 x 20µl
JMJD2 Assay Buffer	1 x 4ml	1 x 8ml
JMJD2 Assay Standard, 50 µg/mL	1 x 10µl	1 x 20µl
JMJD2 Substrate, 50 µg/mL	1 x 60µl	1 x 120µl

## Images



Demonstration of high sensitivity of the KDM4/JMJD2 activity assay achieved by using recombinant KDM4/JMJD2 with ab113462.

**Please note:** All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

## 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