

## Product datasheet

# Histone H3 (K27) Methyltransferase Activity Quantification Assay Kit ab113454

### 画像数 1

#### 製品の概要

製品名	Histone H3 (K27) Methyltransferase Activity Quantification Assay Kit
アッセイタイプ	Quantitative
全工程の試験時間	3h 00m
種交差性	交差種: Mouse, Human 交差が予測される動物種: Mammal

#### 製品の概要

Histone H3 (K27) Methyltransferase Activity Quantification Assay Kit (ab113454) allows the user to measure the activity or inhibition of individual histone methyltransferase that specifically target histone H3 at lysine 27 (*H3K27*). The kit is ready-to-use and provides all the essential components needed to carry out a successful HMT activity/inhibition quantification experiment without the need for radioactivity or any special equipment.

Methylation of histone H3 at lysine 27 (K27) is a mark of gene inactivation and it is thought to mediate heterochromatin formation by forming a binding site for Polycomb. In mammalian cells, the EZ family group of proteins among others are responsible for the addition of one, two or three methyl groups at lysine 27 of histone H3.

#### アプリケーション

適用あり: Functional Studies

#### 製品の特徴

保存方法 Please refer to protocols.

内容	48 tests	96 tests
10X Wash Buffer	1 x 11ml	1 x 22ml
8-Well Assay Strips (with Frame)	6 units	12 units
Adomet	1 x 25µl	1 x 50µl
Biotinylated Substrate, 25 µg/mL	1 x 100µl	1 x 200µl

内容	48 tests	96 tests
Capture Antibody, 100 µg/mL	1 x 25µl	1 x 50µl
Control Enzyme (150 µg/ml)	1 x 10µl	1 x 20µl
Detection Antibody, 100 µg/mL	1 x 10µl	1 x 20µl
Developing Solution	1 x 6ml	1 x 12ml
Histone Assay Buffer	1 x 1.5ml	1 x 3ml
HMT Standard, 10 µg/mL	1 x 10µl	1 x 20µl
Stop Solution	1 x 3ml	1 x 6ml

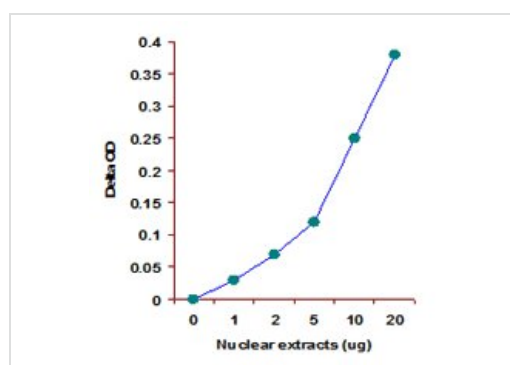
## アプリケーション

Our [Abpromise guarantee](#) covers the use of **ab113454** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

アプリケーション	Abreviews	特記事項
Functional Studies		Use at an assay dependent concentration.

## 画像



Nuclear extracts were prepared from MCF-7 cells using the [ab113474](#) and H3K27 specific histone methyltransferase activity was measured.

Functional Studies - Histone H3 (K27)  
Methyltransferase Activity Quantification Assay Kit  
(ab113454)

**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 <http://www.abcam.co.jp/abpromise> or contact our technical team.

### **Terms and conditions**

---

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