

## Product datasheet

# Histone H3 (tri-methyl K27) Quantification Kit (Fluorometric) ab115073

### 1 References

#### 概述

产品名称	Histone H3 (tri-methyl K27) Quantification试剂盒(Fluorometric)
样品类型	Tissue, Adherent cells, Suspension cells
检测类型	Quantitative
灵敏度	$\geq 0.4$ ng/well
范围	5 ng/well - 2000 ng/well
检测时间	2h 30m
种属反应性	与反应: Mouse, Human 预测可用于: Mammal

#### 产品概述

Methylation of histone H3 at lysine 27 has an important role in transcription repression and it is catalyzed by G9a and Polycomb Groups enzymes such as EZH2. In particular, *H3(tri methyl K27)* is a facultative heterochromatin mark, which promotes the recruitment of polycomb group proteins for gene silencing. Increased global histone H3 (tri methyl K27) is also found to be involved in some pathological processes such as cancer progression.

Histone H3 (tri-methyl K27) Quantification Kit (Fluorometric) (ab115073) allows the user to specifically measure global tri-methylation of histone H3K27 fluorometrically using a variety of mammalian cells including fresh and frozen tissues, cultured adherent and suspension cells.

#### 经测试应用

适用于: Functional Studies

#### 性能

存放说明 Please refer to protocols.

组件	标识符	48 tests	96 tests
10X Wash Buffer		1 x 10ml	1 x 20ml
8-Well Sample Strips (with Frame)		4 units	9 units
8-Well Standard Control Strips	Green Ringed	2 units	3 units
Antibody Buffer		1 x 6ml	1 x 12ml

组件	标识符	48 tests	96 tests
Detection Antibody, 1 mg/mL		1 x 5µl	1 x 10µl
Fluoro-Developer		1 x 12µl	1 x 24µl
Fluoro-Dilution		1 x 4ml	1 x 8ml
Fluoro-Enhancer		1 x 12µl	1 x 24µl
Standard Control (100 µg/mL)		1 x 10µl	1 x 20µl

## 应用

Our [Abpromise guarantee](#) covers the use of **ab115073** in the following tested applications.

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

应用	Ab评论	说明
Functional Studies		Use at an assay dependent concentration.

**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.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