

Product datasheet

Hydroxymethylated DNA Quantification Kit (Fluorometric)  
ab117131

画像数 1

製品の概要

製品名	Hydroxymethylated DNA Quantification Kit (Fluorometric)
サンプルの種類	Saliva, Urine, Plasma, Tissue, Adherent cells, Suspension cells
アッセイタイプ	Quantitative
検出感度	>= 10 pg/well
全工程の試験時間	3h 20m
種交差性	交差種: Mouse, Human

交差が予測される動物種: Plants, Mammal, Fungi ▲

製品の概要

DNA methylation is a covalent modification of the cytosine ring at the 5' position, resulting in 5-*methylcytosine* (5-*mC*). In somatic cells, 5-*mC* is found almost exclusively in the context of paired symmetrical methylation of the dinucleotide CpG. The biological important of 5-*mC* as a major epigenetic modification in phenotype and gene expression has been recognized widely. Quite recently, a novel modified nucleotide, **5-hydroxymethylcytosine (5-hmC)**, has been detected to be abundant in mouse brain and embryonic stem cells. It is a hydroxylated and methylated form of cytosine. Although it is still uncertain its specific role in epigenetics, it is believed it plays a role in DNA demethylation, chromatin remodeling and gene expression regulation.

Hydroxymethylated DNA Quantification Kit (Fluorometric) (ab117131) is a complete set of optimized buffers and reagents to quantify global DNA hydroxymethylation by specifically measuring levels of 5-hydroxymethylcytosine (5-hmC) without cross-reactivity to methylcytosine and unmethylated cytosine.

It can be paired to use with [EpiSeeker methylated DNA Quantification Kit \(Fluorometric\) \(ab117129\)](#), which specifically quantifies **5-mC**, for simultaneously quantification of both methylated and hydroxymethylated DNA.

アプリケーション

適用あり: Functional Studies

製品の特性

保存方法

Please refer to protocols.

内容	48 tests	96 tests
10X Wash Buffer	1 x 14ml	1 x 28ml
8-Well Assay Strips (with Frame)	6 units	12 units
Binding Solution	1 x 5ml	1 x 10ml
Capture Antibody, 1000 µg/mL	1 x 4µl	1 x 8µl
Detection Antibody, 400 µg/mL	1 x 8µl	1 x 16µl
Enhancer Solution	1 x 8µl	1 x 16µl
Fluoro Developer	1 x 8µl	1 x 16µl
Fluoro Dilutor	1 x 4ml	1 x 8ml
Fluoro Enhancer	1 x 8µl	1 x 16µl
Negative Control I, 20 µg/mL	1 x 10µl	1 x 20µl
Negative Control II, 20 µg/mL	1 x 10µl	1 x 20µl
Positive Control, 20 µg/mL	1 x 10µl	1 x 20µl

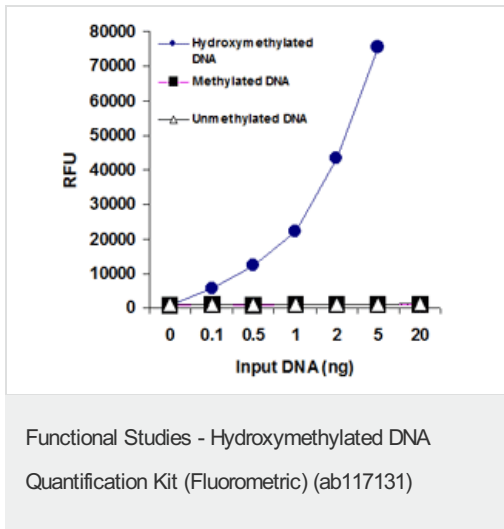
## アプリケーション

Our [Abpromise guarantee](#) covers the use of **ab117131** 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



Demonstration of high sensitivity and specificity of 5-hydroxymethylcytosine detection achieved using ab117131. Synthetic unmethylated DNA (contains only cytosine), methylated DNA (contains only 5-methylcytosine), and hydroxymethylated DNA standard (contains only 5-hydroxymethylcytosine) were added into the assay wells at different concentrations and then measured with ab117131.

**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