

Hydroxymethylated DNA Quantification Kit (Colorimetric)
ab117130

5 References 画像数 4

製品の概要

製品名	Hydroxymethylated DNA Quantification Kit (Colorimetric)
検出方法	Colorimetric
サンプルの種類	Saliva, Urine, Serum, Plasma, Tissue, Adherent cells, Suspension cells
アッセイタイプ	Quantitative
検出感度	>= 40 pg/well
全工程の試験時間	3h 45m
製品の概要	<p>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.</p> <p>Abcam's Hydroxymethylated DNA Quantification Kit (Colorimetric) (ab117130) 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.</p> <p>It can be paired to use with Methylated DNA Quantification Kit (Colorimetric) (ab117128), which specifically quantifies 5-mC, for simultaneously quantification of both methylated and hydroxymethylated DNA.</p>
試験プラットフォーム	Microplate reader

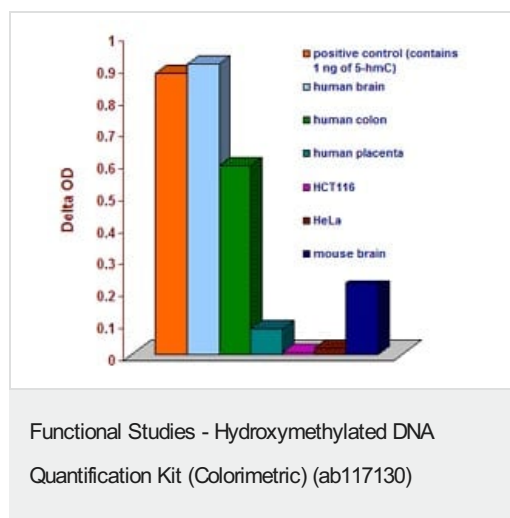
製品の特性

保存方法 Please refer to protocols.

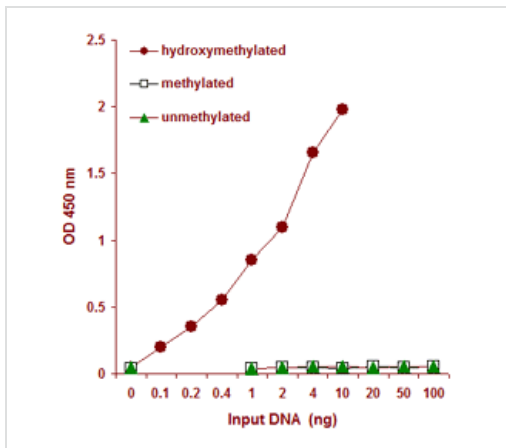
内容	48 tests	96 tests
10X Wash Buffer	1 x 14ml	1 x 28ml

内容	48 tests	96 tests
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
Developer Solution	1 x 5ml	1 x 10ml
Enhancer Solution	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
Stop Solution	1 x 5ml	1 x 10ml

画像

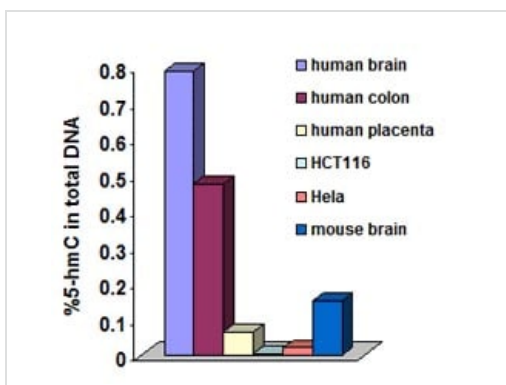


Demonstration of a high abundance of 5-hydroxymethylcytosine in human brain and colon tissues. 200 ng of DNA isolated from different tissues or cells were added into the assay wells and the hydroxymethylated fraction of DNA was measured using the ab117130. OD intensity measured in variety of DNA samples.



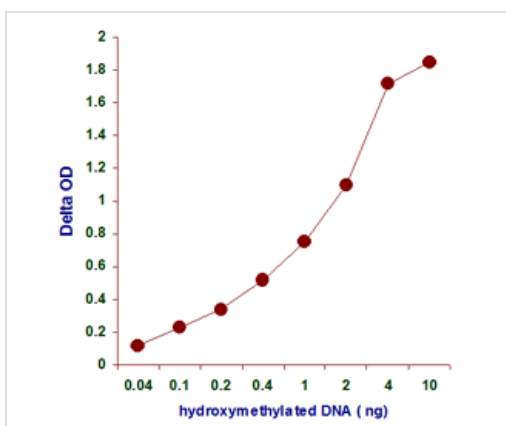
Functional Studies - Hydroxymethylated DNA
Quantification Kit (Colorimetric) (ab117130)

Demonstration of high specificity of 5-hydroxymethylcytosine detection achieved using ab117130. 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 ab117130.



Functional Studies - Hydroxymethylated DNA
Quantification Kit (Colorimetric) (ab117130)

Demonstration of a high abundance of 5-hydroxymethylcytosine in human brain and colon tissues. 200 ng of DNA isolated from different tissues or cells were added into the assay wells and the hydroxymethylated fraction of DNA was measured using ab117130. Percentage of hydroxymethylated DNA in various tissues and cells.



Functional Studies - Hydroxymethylated DNA
Quantification Kit (Colorimetric) (ab117130)

Demonstration of high sensitivity of 5-hydroxymethylcytosine detection achieved using ab117130. Synthetic hydroxymethylated DNA was added into the assay wells at different concentrations and then measured with the ab117130.

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.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