

Methylated DNA Quantification Kit (Colorimetric) ab117128

30 References 画像数 1

製品の概要

製品名	Methylated 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	>= 0.2 ng/well
全工程の試験時間	4h 00m
種交差性	交差種: Species independent
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. Quantification of 5-mC content or global methylation status in cells such as tumor cells could provide very useful information for detection and analysis of cancer.</p> <p>Methylated DNA Quantification Kit (Colorimetric) (ab117128) is a complete set of optimized buffers and reagents to quantify global DNA methylation by specifically measuring levels of 5-methylcytosine (5-mC) in a microplate-based format.</p> <p>It can be paired to use with <u>Abcam's Hydroxymethylated DNA Quantification Kit (Colorimetric) (ab117130)</u>, which specifically quantifies 5-hmC, for simultaneous quantification of both methylated and hydroxymethylated DNA.</p>
試験プラットフォーム	Microplate reader

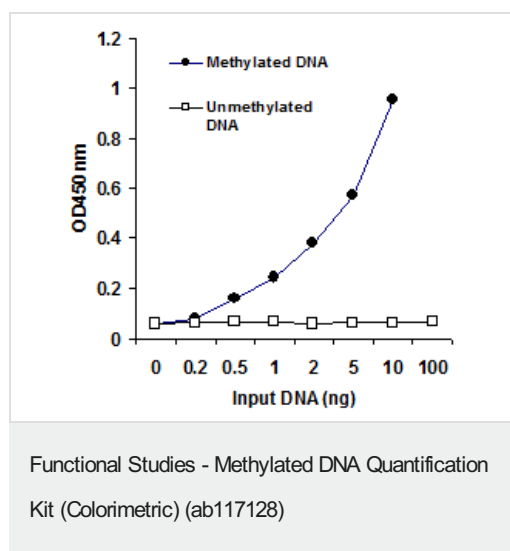
製品の特性

保存方法 Please refer to protocols.

内容	48 tests	96 tests
10X Wash Buffer	1 x 14ml	1 x 28ml
8-Well Assay Strips (with Frame)	6 units	12 units
Binding Solution	1 x 5ml	1 x 10ml

内容	48 tests	96 tests
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, 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 high sensitivity and specificity of methylated DNA detection achieved using ab117128. Synthetic unmethylated DNA (contains 50% of cytosine) and methylated DNA (contains 50% of 5-methylcytosine) were added into the assay wells at different concentrations and then measured with ab117128.

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