

Product datasheet

Methylated DNA Quantification Kit (Fluorometric) ab117129

★★★★☆ [1 Abreviews](#) [2 References](#) [1 Image](#)

Overview

Product name	Methylated DNA Quantification Kit (Fluorometric)
Detection method	Fluorescent
Sample type	Saliva, Urine, Plasma, Tissue, Adherent cells, Suspension cells
Assay type	Quantitative
Sensitivity	>= 50 pg/well
Assay time	4h 00m
Species reactivity	Reacts with: Species independent
Product overview	<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 importance 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>Abcam's Methylated DNA Quantification Kit (Fluorometric) (ab117129) 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 Hydroxymethylated DNA Quantification Kit (Fluorometric) (ab117131), which specifically quantifies 5-hmC, for simultaneous quantification of both methylated and hydroxymethylated DNA.</p>
Platform	Microplate reader

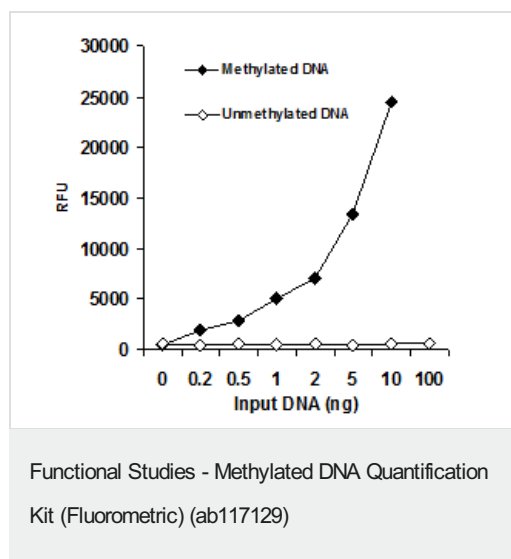
Properties

Storage instructions Please refer to protocols.

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

Components	48 tests	96 tests
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, 20 µg/mL	1 x 10µl	1 x 20µl
Positive Control, 20 µg/mL	1 x 10µl	1 x 20µl

Images



Demonstration of high sensitivity and specificity of methylated DNA detection achieved using ab117129. 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 ab117129.

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.com/abpromise> or contact our technical team.

Terms and conditions

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