


DNMT3B Assay Kit ab113471

3 References 1 图像

概述

产品名称	DNMT3B Assay试剂盒
检测时间	3h 00m
种属反应性	与反应: Mouse, Human 预测可用于: Mammals 
产品概述	<p>DNMT3B (DNA (cytosine-5)-methyltransferase 3B) belongs to the DNMT family, it is required for genome <i>de novo</i> methylation and it is essential for establishment of DNA methylation patterns during development.</p> <p>DNMT3B Assay Kit allows the user to quickly and consistently measure the amount of DNMT3B. The kit is ready-to-use and provides all the essential components needed to carry out a successful DNMT3B assay experiment. DNMT3B Assay Kit is suitable for quantitative measurement of DNMT3B amounts from fresh tissue and cultured cells of human and mouse.</p>
平台	Microplate reader

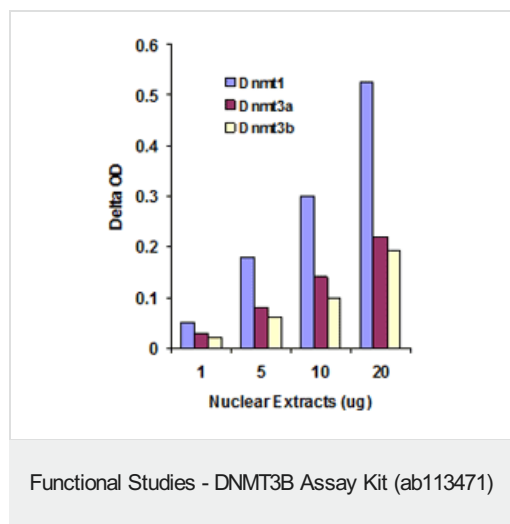
性能

存放说明 Please refer to protocols.

组件	48 tests	96 tests
10X Wash Buffer	1 x 12ml	1 x 25ml
8-Well Assay Strips (with Frame)	6 units	12 units
Assay Buffer	1 x 5ml	1 x 10ml
Blocking Buffer	1 x 10ml	1 x 20ml
Capture Antibody, 100 µg/mL	1 x 8µl	1 x 16µl
Detection Antibody, 200 µg/mL	1 x 10µl	1 x 20µl
Developing Solution	1 x 6ml	1 x 12ml

组件	48 tests	96 tests
DNMT3B Standard, 20 µg/mL	1 x 16µl	1 x 30µl
Stop Solution	1 x 6ml	1 x 11ml

图片



Nuclear extracts were prepared from MCF-7 cells using the Nuclear Extraction Kit ([ab113474](#)) and DNMT3B activity was measured using ab113471. DNMT1 and DNMT3A activities were measured using [ab113469](#) DNMT1 Assay Kit and [ab113470](#) DNMT3A Assay Kit.

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.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