

RNA Bisulfite Conversion Kit ab185911

概述

产品名称	RNA Bisulfite Conversion试剂盒
灵敏度	5 ng
检测时间	3h 00m
实验步骤	Multiple steps standard assay
产品概述	The RNA Bisulfite Conversion Kit (ab185911) is a complete set of optimized reagents required for fast bisulfite conversion on a RNA sample. The converted RNA obtained with the RNA Bisulfite Conversion Kit is suitable for various downstream RNA methylation analysis including methylation specific RT-PCR, MS-HRM, and bisulfite RNA-sequencing (e.g., pyrosequencing and deep-sequencing).

说明	<p>DNA cytosine methylation occurs by the covalent addition of a methyl group at the 5-carbon of the cytosine ring by DNA methyltransferases, resulting in 5-methylcytosine (5-mC). This process has been well studied and is generally associated with repression of gene expression. It was also observed that in humans, 5-mC occurs in various RNA molecules including tRNAs, rRNAs, mRNAs and non-coding RNAs (ncRNAs). At least 10,275 5-mC candidate sites were discovered in mRNAs and ncRNAs, which covered 10.6% of the total cytosine residues in the transcriptome. 5-mC seems to be enriched in some classes of ncRNA, but relatively depleted in mRNAs. However, the majority (83%) of their candidate sites were found in mRNAs. Within these transcripts 5-mC appears to be depleted within protein coding sequences but enriched in 5' and 3' UTRs. Two different methyltransferases, NSUN2 and Dnmt2 are known to catalyze the 5-mC modification in eukaryotic RNAs. Recent data strongly suggest that RNA cytosine methylation affects the regulation of various biological processes such as RNA stability and mRNA translation. Furthermore, loss of 5-mC in vault RNAs causes aberrant processing into Argonaute-associated small RNA fragments that can function as microRNAs. Thus, impaired processing of vault ncRNA may contribute to the etiology of human disorders related to NSUN2-deficiency.</p>
----	---

性能

存放说明	Please refer to protocols.
------	----------------------------

组件	50 tests
Control Primer F	1 x 10µl
Control Primer R	1 x 10µl

组件	50 tests
Conversion Buffer	1 x 8ml
Conversion Powder	1 x 5 vials
Desulphonation Solution	1 x 300µl
Elution Buffer	1 x 1.2ml
F-Collection Tube	1 x 50 units
F-Spin Column	1 x 50 units
NA Binding Solution	1 x 13ml

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