

MBD2 Binding Activity Quantification Assay Kit ab113473

製品の概要

製品名	MBD2 Binding Activity Quantification Assay Kit
アッセイタイプ	Quantitative
全工程の試験時間	3h 00m
種交差性	交差種: Mouse, Human 交差が予測される動物種: Mammals

製品の概要

MBD2 (methyl-CpG binding domain protein 2) belongs to a family of MBD domain containing proteins. This family of proteins associates with heterochromatin by their interaction with methylated DNA at CpG islands, and appear to 'read' and 'translate' the DNA methylation signal into transcriptional repression by recruiting silencing complexes and histone deacetylases among other proteins. MBD2 is the methyl-binding component of the MeCP1 complex, which allows MBD2 to target HDAC/chromatin remodeling activity to methylated templates.

MBD2 Binding Activity Quantification Assay Kit (ab113473) allows the user to measure binding activity of MBD2 to methylated DNA in human cells/tissues. It can also be used to determine inhibitors of MBD2 activity.

製品の特性

保存方法 Please refer to protocols.

内容	48 tests	96 tests
10X Wash Buffer	1 x 12ml	1 x 25ml
8-Well Assay Strips (with Frame)	6 units	12 units
Affinity Antibody, 100 µg/mL	1 x 25µl	1 x 50µl
Assay Buffer	1 x 1.8ml	1 x 4ml
Detection Antibody, 200 µg/mL	1 x 10µl	1 x 20µl
Developing Solution	1 x 6ml	1 x 12ml
Methylated DNA, 25 µg/mL	1 x 50µl	1 x 100µl

内容	48 tests	96 tests
Stop Solution	1 x 3ml	1 x 6ml

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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