

ChIP Kit (FFPE samples) ab270815

製品の概要

製品名 ChIP Kit (FFPE samples)

製品の概要 ChIP Kit (FFPE samples) (ab270815) provides a unique and robust ChIP protocol suitable for investigation of proteins associated with chromatin from FFPE samples and provides high quality yields of DNA. This solution makes traditionally difficult work with FFPE samples simpler and safer. The protocol involves easy and quick and non-toxic deparaffinization, followed by chromatin extraction and shearing by sonication. The kit contains everything you need for start-to-finish chromatin immunoprecipitation with an optimized protocol to make ChIP from FFPE samples simple yet reliable. In addition, the kit has been thoroughly validated and optimized for histones and transcription factors.

This kit contains enough reagents to perform 10 individual chromatin preparation from FFPE sections, up to 24 immunoprecipitations and 50 DNA purifications (from 10 input, 24 IPs and 10 shearing assessment samples).

特記事項

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances. It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

製品の特性

保存方法 Please refer to protocols.

内容	24 tests
5% BSA (DNA free)	1 x 144µl
5x ChIP Buffer C1b	1 x 6.8ml
ChIP DNA Binding Buffer	1 x 50ml
ChIP-seq grade water	1 x 26.6ml
DNA Elution Buffer	1 x 10ml
DNA Wash Buffer	1 x 6ml

内容	24 tests
Elution Buffer E1	1 x 7.8ml
Elution Buffer E2	1 x 320µl
Filters	10 units
Lysis Buffer L1b	1 x 6ml
Lysis Buffer L2	1 x 6ml
Protease inhibitor cocktail	1 x 80µl
Protein A-coated magnetic beads	1 x 480µl
Shearing Buffer S1b	1 x 6ml
Spin columns	50 units
Wash buffer W1	1 x 8.4ml
Wash buffer W2	1 x 8.4ml
Wash buffer W3	1 x 8.4ml
Wash Buffer W4	1 x 8.4ml

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