

ChIP Kit (Magnetic, qPCR) ab270816

1 References

製品の概要

製品名	ChIP Kit (Magnetic, qPCR)
製品の概要	ChIP Kit (Magnetic, qPCR) (ab270816) is a highly optimized and validated solution for ChIP-qPCR assays for either histones or transcription factors. The complete kit contains everything you need for start-to-finish ChIP including all validated buffers and reagents for chromatin shearing, immunoprecipitation, and DNA isolation for exceptional ChIP-qPCR results.

This kit uses a unique and fast method for DNA isolation and decrosslinking within 30 minutes compared to the standard 4 hours. Overall, easy to use and rapid protocol allows to receive the results within 20 hours with 4 hours hands-on time.

This kit contains enough reagents to perform 4 chromatin preparations, 24 chromatin immunoprecipitations and 32 DNA isolations.

Sample size:

Histones -

Cells (100,000 cells to 1,000,000 cells per IP)

Tissues (1.5 mg to 5 mg of tissue per IP)

Transcription factors -

Cells (4,000,000 cells per IP)

Tissues (30 mg of tissue per IP)

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 784µl
5x ChIP Buffer C1b	1 x 6.9ml
ChIP-seq grade water	1 x 26.6ml
Protein A-coated magnetic beads	1 x 720µl
DNA isolation buffer (DIB)	1 x 3.2ml
Elution Buffer E1	1 x 220µl
Elution Buffer E2	1 x 20µl
Fixation buffer	1 x 8ml
Glycine	1 x 8.8ml
Lysis Buffer L1b	1 x 100ml
Lysis Buffer L2	1 x 60ml
Protease inhibitor cocktail	1 x 80µl
Proteinase K	1 x 32µl
Rabbit IgG	1 x 8µg
Shearing Buffer S1b	1 x 9ml
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