



Product datasheet

ChIP Kit (Histones, ChIP-seq) ab270812

概述

产品名称	ChIP试剂盒(Histones, ChIP-seq)
产品概述	<p>ChIP Kit (Histones, ChIP-seq) (ab270812) provides a robust ChIP protocol suitable for the investigation of histone modifications within chromatin from cells and tissues. The protocol involves protein-DNA cross-linking with formaldehyde, followed by cell lysis and fragmentation of the cross-linked chromatin. The subsequent immunoprecipitation of chromatin is performed with an antibody (user supplied) specific to a target histone modification. Protein A magnetic beads are used to isolate the protein-DNA complexes of interest. The immunoprecipitated DNA is eluted and purified using unique IPure magnetic beads. The eluted DNA is ideal for ChIP-seq library preparation. Additionally, the kit includes a negative (IgG) and a positive (H3K4me3) control antibody and primer pairs amplifying H3K4me3 positive and negative loci.</p> <p><b>Sample size:</b></p> <p>Cells (100,000 cells to 1,000,000 cells per IP)</p> <p>Tissues (1.5 mg to 5 mg of tissue per IP)</p>

说明	<p>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.</p> <p>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.</p>
----	--

性能

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

组件	24 tests	10 tests
5% BSA (DNA free)	1 x 144µl	1 x 60µl
5x ChIP Buffer C1	1 x 2.2ml	1 x 950µl
Buffer C	1 x 1.6ml	1 x 700µl
Carrier	1 x 64µl	1 x 28µl

组件	24 tests	10 tests
ChIP-seq grade antibody H3K4me3	1 x 8µg	1 x 8µg
ChIP-seq grade GAPDH TSS primer pair	1 x 96µl	1 x 42µl
ChIP-seq grade Myoglobin exon 2 primer pair	1 x 96µl	1 x 42µl
ChIP-seq grade water	1 x 7ml	1 x 3ml
Elution Buffer E1	1 x 3.4ml	1 x 1.5ml
Elution Buffer E2	1 x 144µl	1 x 64µl
Glycine	1 x 400µl	1 x 200µl
Lysis Buffer L1	1 x 40ml	1 x 20ml
Lysis Buffer L2	1 x 40ml	1 x 20ml
Protease inhibitor cocktail	1 x 76µl	1 x 35µl
Protein A-coated magnetic beads	1 x 480µl	1 x 200µl
Purification Beads	1 x 320µl	1 x 140µl
Rabbit IgG	1 x 8µg	1 x 8µg
Shearing Buffer S1	1 x 7.2ml	1 x 3.6ml
Wash Buffer 1 w/o iso-propanol	1 x 2ml	1 x 700µl
Wash Buffer 2 w/o iso-propanol	1 x 2ml	1 x 700µl
Wash Buffer W1	1 x 8.4ml	1 x 3.5ml
Wash Buffer W2	1 x 8.4ml	1 x 3.5ml
Wash Buffer W3	1 x 8.4ml	1 x 3.5ml
Wash Buffer W4	1 x 8.4ml	1 x 3.5ml

**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