

ChIP Kit (Transcription factors, ChIP-seq) ab270813

4 图像

概述

产品名称

ChIP试剂盒(Transcription factors, ChIP-seq)

产品概述

ChIP Kit (Transcription factors, ChIP-seq) (ab270813) is a highly validated solution for robust transcription factor and other non-histone proteins ChIP-seq results and contains everything you need for start-to-finish ChIP prior to Next-Generation Sequencing. This complete solution contains all buffers and reagents for cell lysis, chromatin shearing, immunoprecipitation, and DNA purification. This kit contains positive and negative control antibodies (CTCF and IgG, respectively) as well as positive and negative control PCR primers pairs (H19 and Myoglobin exon 2, respectively).

Sample size:

Cells (4,000,000 cells per IP)

Tissues (30 mg of tissue per IP)

说明

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.

性能

存放说明

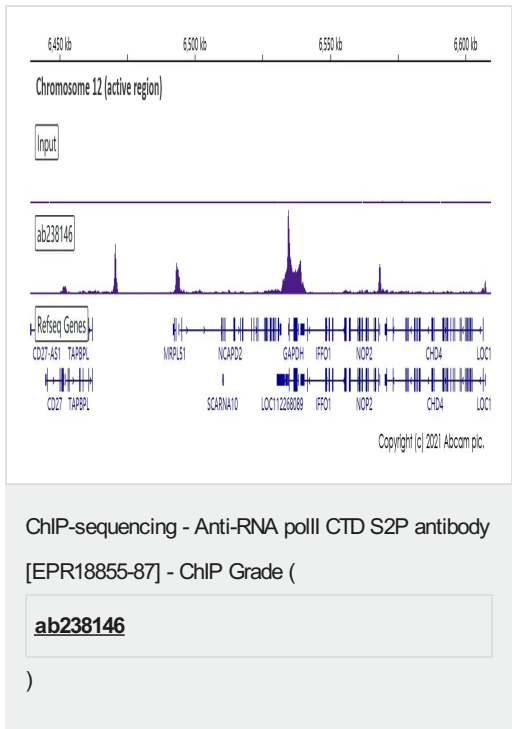
Please refer to protocols.

组件	24 tests	10 tests
5% BSA (DNA free)	1 x 800µl	1 x 380µl
5x ChIP Buffer C1b	1 x 6.9ml	1 x 3.4ml
Buffer C	1 x 1.6ml	1 x 700µl
Carrier	1 x 72µl	1 x 32µl

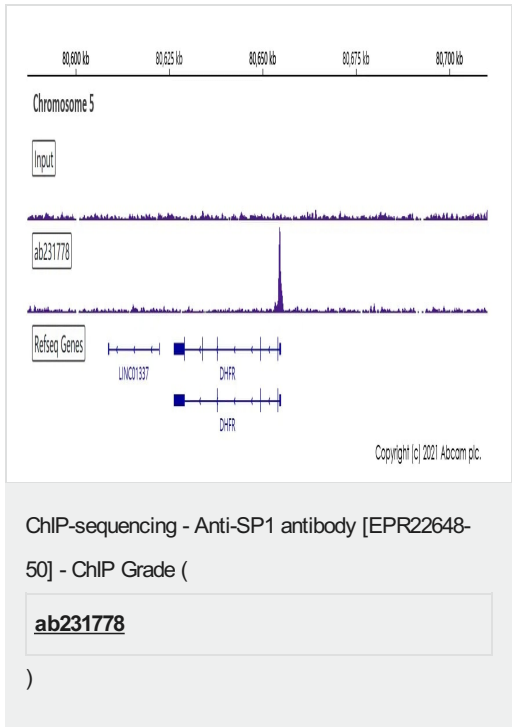
组 件	24 tests	10 tests
ChIP-seq grade CTCF antibody	1 x 8μg	1 x 4μg
ChIP-seq grade H19 imprinting control region primer pair (human)	1 x 96μl	1 x 42μl
ChIP-seq grade Myoglobin exon 2 primer pair (human)	1 x 96μl	1 x 42μl
ChIP-seq grade water	1 x 26.6ml	1 x 14ml
Elution Buffer E1	1 x 3.4ml	1 x 1.5ml
Elution-buffer E2	1 x 144μl	1 x 64μl
Fixation buffer	1 x 8ml	1 x 4ml
Glycine	1 x 8.8ml	1 x 4.4ml
Lysis Buffer L1b	1 x 100ml	1 x 50ml
Lysis Buffer L2	1 x 60ml	1 x 30ml
Protease inhibitor cocktail	1 x 80μl	1 x 38μl
Protein A-coated magnetic beads	1 x 720μl	1 x 300μl
Purification Beads	1 x 400μl	1 x 180μl
Rabbit IgG	1 x 8μg	1 x 4μg
Shearing Buffer S1b	1 x 6.7ml	1 x 3.4ml
Wash buffer 1 w/o iso-propanol	1 x 2ml	1 x 900μl
Wash buffer 2 w/o iso-propanol	1 x 2ml	1 x 900μl
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

## 细胞定位

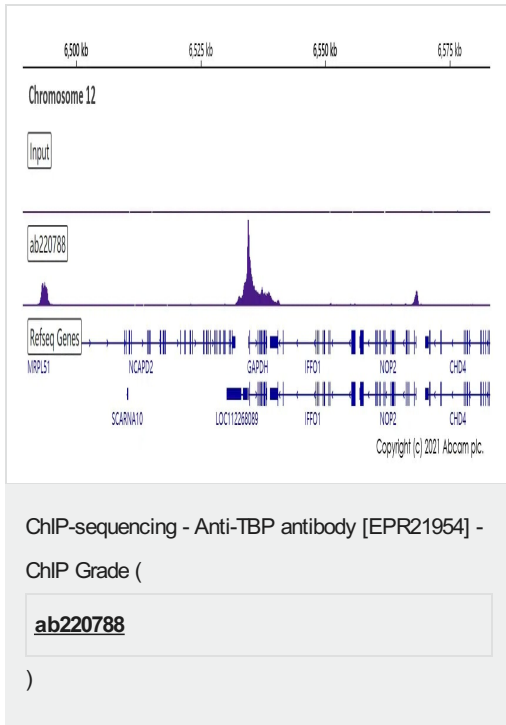
RNA polymerase II CTD repeat YSPTSPS: Nucleus. TATA binding protein TBP: Nucleus. SP1: Nucleus. Cytoplasm. Nuclear location is governed by glycosylated/phosphorylated states. Insulin promotes nuclear location, while glucagon favors cytoplasmic location. CTCF: Nucleus > nucleoplasm. Chromosome. Chromosome > centromere. May translocate to the nucleolus upon cell differentiation. Associates with both centromeres and chromosomal arms during metaphase. Associates with the H19 ICR in mitotic chromosomes. May be preferentially excluded from heterochromatin during interphase.



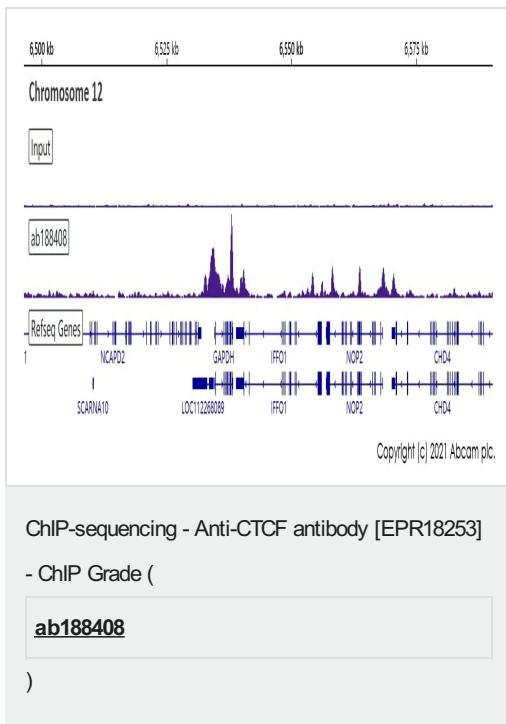
Chromatin was prepared from HeLa cells. Cells were fixed with 1% formaldehyde for 10 minutes. ChIP was performed with 30 µg of chromatin and 4 µg of **ab238146**. ChIP DNA was sequenced on the Illumina NextSeq 500 to a depth of 30 million reads. ChIP-Seq validation performed with ChIP-Kit Transcription Factors ChIP-Seq (ab270813).



Chromatin was prepared from HeLa cells. Cells were fixed with 1% formaldehyde for 10 minutes. ChIP was performed with 30 µg of chromatin and 4 µg of **ab231778**. ChIP DNA was sequenced on the Illumina NextSeq 500 to a depth of 30 million reads. ChIP-Seq validation performed with ChIP-Kit Transcription Factors ChIP-Seq (ab270813).



Chromatin was prepared from HeLa cells. Cells were fixed with 1% formaldehyde for 10 minutes. ChIP was performed with 30 µg of chromatin and 4 µg of **ab220788**. ChIP DNA was sequenced on the Illumina NextSeq 500 to a depth of 30 million reads. ChIP-Seq validation performed with ChIP-Kit Transcription Factors ChIP-Seq (ab270813).



Chromatin was prepared from HeLa cells. Cells were fixed with 1% formaldehyde for 10 minutes. ChIP was performed with 30 µg of chromatin and 4 µg of **ab188408**. ChIP DNA was sequenced on the Illumina NextSeq 500 to a depth of 30 million reads. ChIP-Seq validation performed with ChIP-Kit Transcription Factors ChIP-Seq (ab270813).

### **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