

ChIP Kit - Plants ab117137

8 References 2 图像

概述

产品名称	ChIP试剂盒- Plants
检测类型	Quantitative
检测时间	6h 00m
种属反应性	与反应: Plants
产品概述	<p>Protein-DNA interactions play a critical role for cellular functions such as signal transduction, gene transcription and epigenetic silencing. In plants, interactions between the DNA-binding proteins and cognate promoter sequences are primary determinants in establishing spatial and temporal expression patterns of gene that affect homeostasis, development, and adaptation.</p> <p>Abcam's ChIP Kit - Plants (ab117137) offers an advantageous tool for identifying direct genome-wide associations between specific regulatory proteins and their target genes in plant cells within 6 hours. The kit is suitable for combining the specificity of immunoprecipitation with qualitative and quantitative PCR, MS-PCR, DNA sequencing and southern blot, as well as DNA microarray.</p> <p>This kit is designed for 24 or 48 ChIP reactions, not for 24 or 48 samples. The standard protocol of the kit allows for performing 8 reactions with one sample. For testing more samples, the amount of each sample should be reduced. The amount of each reagent used for chromatin preparation should be also proportionally reduced.</p>

**说明**

**ChIP assay products and guides**

Find more [ChIP assay / chromatin immunoprecipitation](#) resources and products, [ChIP antibody](#) products, and other [ChIP assay kits and related reagents](#).

**经测试应用**

**适用于:** ChIP

性能

**存放说明**

Please refer to protocols.

组件	24 tests	48 tests
100X Protease Inhibitor Cocktail	1 x 25µl	1 x 50µl
5X Lysis Buffer I	1 x 12ml	1 x 24ml

组件	24 tests	48 tests
8-Well Assay Strips (with Frame)	3 units	6 units
8-Well Strip Caps	3 units	6 units
Antibody Buffer	1 x 15ml	1 x 30ml
Anti-H3K9me2 (1 mg/mL)	1 x 5µl	1 x 8µl
Binding Buffer	1 x 5ml	1 x 8ml
ChIP Dilution Buffer	1 x 2ml	1 x 6ml
DNA Release Buffer	1 x 2ml	2 x 2ml
Elution Buffer	1 x 0.6ml	1 x 1.2ml
F-Collection Tube	30 units	50 units
F-Spin Column	30 units	50 units
Lysis Buffer II	1 x 3ml	1 x 6ml
Lysis Buffer III	1 x 2ml	1 x 4ml
Lysis Buffer IV	1 x 1.5ml	1 x 5ml
Normal Mouse IgG (1 mg/mL)	1 x 10µl	1 x 10µl
Proteinase K (10 mg/mL)	1 x 25µl	1 x 50µl
Reverse Buffer	1 x 2ml	2 x 2ml
Wash Buffer	1 x 28ml	2 x 28ml

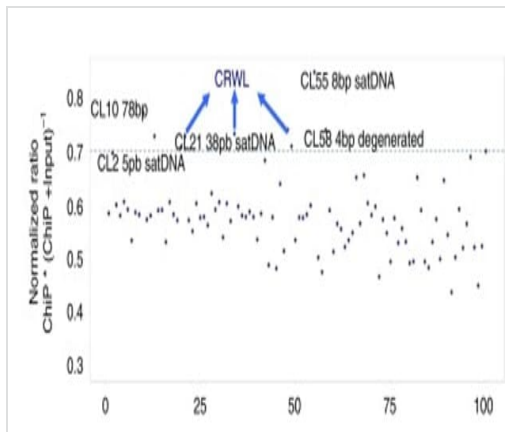
## 应用

**The Abpromise guarantee**      **Abpromise™** 承诺保证使用 ab117137 于以下的经测试应用

“应用说明”部分 下显示的仅为推荐的起始稀释度；实际最佳的稀释度/浓度应由使用者检定。

应用	Ab 评论	说明
ChIP		Use at an assay dependent concentration.

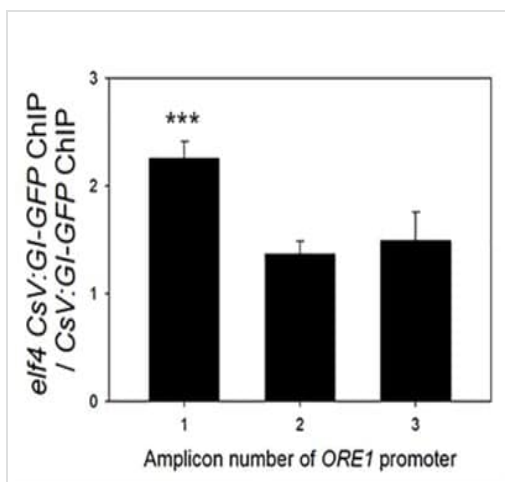
## 图片



ChIP - ChIP Kit - Plants (ab117137)

Image from Hufnagel et al., Nat Commun.;11(1):492. doi: 10.1038/s41467-019-14197-9. Reproduced under the Creative Commons license <https://creativecommons.org/licenses/by/4.0/>

Repeated elements abundance in White Lupin genome. Chromatin immunoprecipitation experiments were done with ChIP Kit - Plants (ab117137). Sonicated chromatin-DNA ranging from 200-1000 bp was immunoprecipitated using anti-LalbCENH3. LalbCENH3-ChIPseq reads mapped against the first 100 RepeatExplorer clusters of the White Lupin genome. The main centromeric sequences found in LalbCENH3-ChIPseq are highlighted.



ChIP - ChIP Kit - Plants (ab117137)

Image from Kim et al., Front Plant Sci.;11:589707. doi: 10.3389/fpls.2020.589707. Reproduced under the Creative Commons license <https://creativecommons.org/licenses/by/4.0/>

ORE1 promoter-binding activity of GIGANTEA in the elf4 mutant relative to that in the wild type at amplicons 1, 2, and 3. ChIP assays were performed using the ChIP Kit-Plants (ab117137). Chromatin was immunoprecipitated using anti-GFP antibody ([ab290](#))-bound assay plate for 90 min. The resulting immunoprecipitated DNA was subjected to qPCR to examine the enrichment of target genes. Three biological replicates were performed. Asterisks indicate significant differences (\* $p < 0.05$ , \*\* $p < 0.01$ , \*\*\* $p < 0.005$ ; Student's t-test).

**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