

15 References 画像数 2

製品の概要

製品名	ChIP Kit - One Step
アッセイタイプ	Quantitative
全工程の試験時間	5h 00m
製品の概要	Abcam's ChIP Kit - One Step (ab117138) enables the user to enrich a chromatin fraction containing specific DNA sequences from isolated chromatin to be able to investigate the interaction between proteins and DNA. For best result, we recommend the use of Chromatin Extraction Kit (ab117152) for sample preparation.

Protein-DNA interactions play a critical role for cellular functions such as signal transduction, gene transcription and epigenetic silencing. In mammalian cells, 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.

特記事項

Primers

The GAPDH primers provided with the kit are for the human sequence. If using the kit with a difference species, GAPDH primers for that species will need to be acquired.

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**.

アプリケーション

適用あり: ChIP

製品の特長

保存方法

Store at +4°C. Please refer to protocols.

内容	ラベル	48 tests	96 tests
10X Wash Buffer		1 x 10ml	1 x 20ml
8-Well Assay Strips (with Frame)		6 units	12 units
8-Well Strip Caps		6 units	12 units

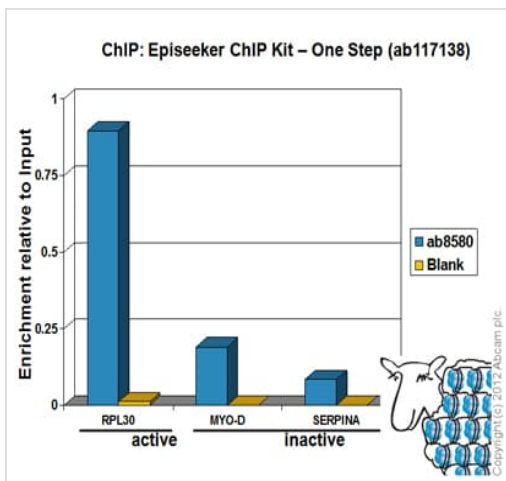
内容	ラベル	48 tests	96 tests
96-Well PCR Plate		1 unit	1 unit
Adhesive Covering Film		1 unit	2 units
Anti-RNA Polymerase II (1 mg/ml)		1 x 8µl	1 x 16µl
ChIP Buffer		1 x 6ml	1 x 12ml
DNA Release Buffer	ECP3	1 x 14ml	1 x 28ml
GAPDH Primer - Forward (20 µM)		1 x 8µl	1 x 16µl
GAPDH Primer - Reverse (20 µM)		1 x 8µl	1 x 16µl
Non-Immune IgG (1 mg/ml)		1 x 10µl	1 x 20µl
Proteinase K (10 mg/mL)		1 x 55µl	1 x 110µl
Sonication Frame		1 unit	1 unit

アプリケーション

The Abpromise guarantee **Abpromise保証は、** 次のテスト済みアプリケーションにおけるab117138の使用に適用されます
 アプリケーションノートには、推奨の開始希釈率がありますが、適切な希釈率につきましてはご検討ください。

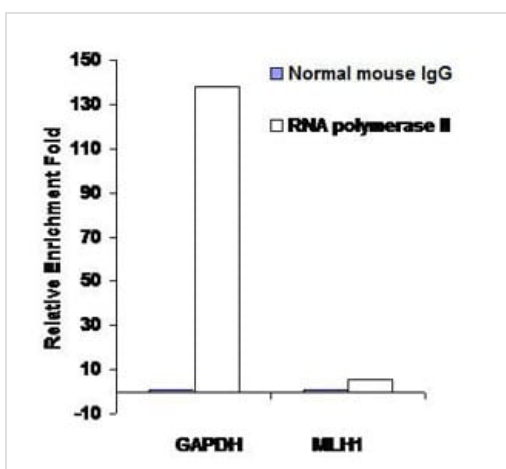
アプリケーション	Abreviews	特記事項
ChIP		Use at an assay dependent concentration. The GAPDH primers provided with the kit are for the human sequence. If using the kit with a different species, GAPDH primers for that species will need to be acquired.

画像



ChIP - ChIP Kit - One Step (ab117138)

Chromatin was prepared from HeLa cells following **Abcam's X-ChIP protocol**. Cells were fixed with formaldehyde for 10 minutes. ChIP was performed with **ab8580 (Histone H3 tri-methyl K4)** using ab117138. No antibody was added to the blank (yellow). The immunoprecipitated DNA was quantified by real time PCR (Taqman approach). Primes and probes are located in the first Kb of the transcribed region.



ChIP - ChIP Kit - One Step (ab117138)

The analysis of enrichment of RNA polymerase II in GAPDH and MLH1 promoters by Abcam's ChIP Kit - One Step, with chromatin extract prepared from formaldehyde fixed colon cancer cells. Captured DNA was used for analyzing levels of RNA polymerase II enriched in the GAPDH and MLH1 promoters.

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