

ChIP Kit ab500

★★★★★ 2 Abreviews 36 References 画像数 4

製品の概要

製品名 ChIP Kit

サンプルの種類 Adherent cells, Suspension cells

製品の概要 ChIP kit ab500 provides a protocol and reagents for running ChIP assays including:

- cell lysis and chromatin extraction
- chromatin shearing and DNA fragment length analysis
- immunoprecipitation and DNA purification

DNA produced using the kit can be analyzed using qPCR.

The kit has been validated for ChIP assays with mammalian samples.

特記事項

This kit uses Protein A sepharose beads for antibody pulldown. See table below for Protein A and Protein G binding affinities with antibodies from commonly used species. For other species of antibody, consult the table in the protocol booklet.

Species raised in	Isotype	Protein A binding affinity	Protein G binding affinity
Rabbit	All isotypes	+++	++
Goat	All isotypes	-	++
	IgG1	+	+++
	IgG2a	+++	+++
	IgG2b	++	++
Mouse	IgG3	+	+
	IgM	Use anti-mouse IgM	

アプリケーション

適用あり: ChIP
適用なし: CHIPseq

製品の特性

保存方法

Please refer to protocols.

バッファー

Kit Constituents: Please see protocol.

内容	1 kit
Buffer A	1 x 10ml
Buffer B	1 x 30ml
1.25M Glycine	1 x 10ml
Buffer C	1 x 30ml
Buffer D	1 x 3ml
ChIP Buffer 5X	1 x 84ml
DNA Purifying Slurry	1 x 3ml
ab1791 - H3 antibody	1 x 25µg
PCR-Grade Water	1 x 10.2ml
Protease Inhibitors	1 tablet
Protein A Beads Unblocked	1 x 960µl
Proteinase K	1 x 30µl

アプリケーション

Our [Abpromise guarantee](#) covers the use of **ab500** in the following tested applications.

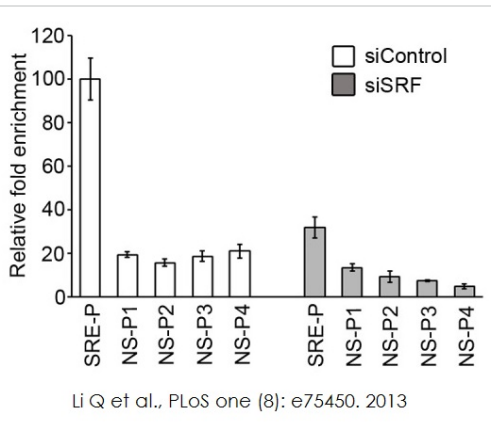
The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

アプリケーション	Abreviews	特記事項
ChIP		Use at an assay dependent concentration.

追加情報

Is unsuitable for CHIPseq.

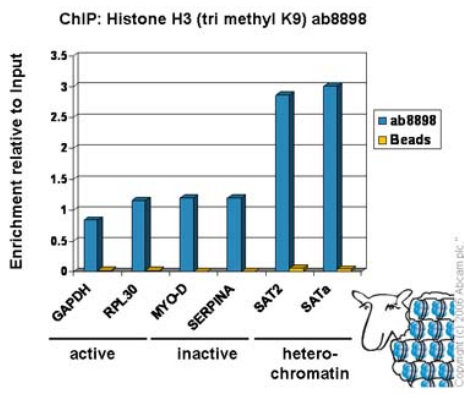
画像



ChIP

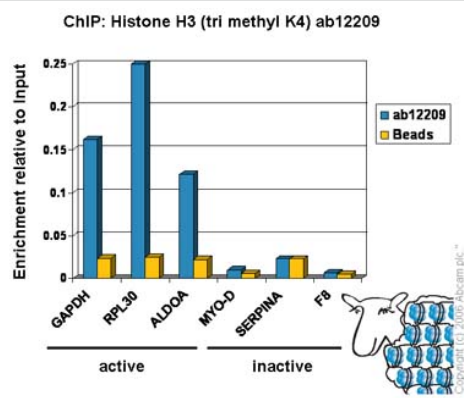
Li Q., PLoS One 8(9). Fig 2c. doi: 10.1371/journal.pone.0075470.

Chromatin immunoprecipitation assay was performed to define the interaction of serum response factor (SRF) with the intragenic serum response element (SRE) regulatory motif in mouse using ab500 ChIP Kit



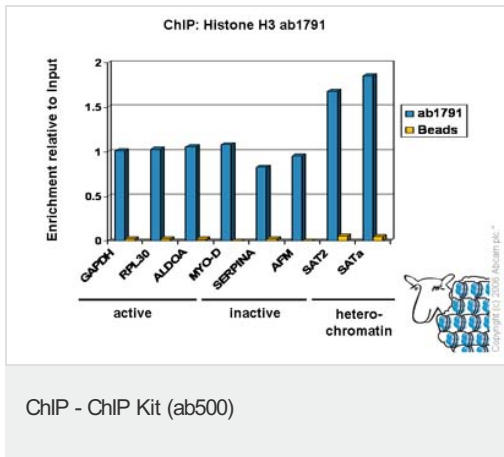
ChIP - ChIP Kit (ab500)

Chromatin immunoprecipitation using ab500 ChIP Kit and Histone H3 (tri methyl K9) antibody (ab8898). Chromatin was prepared from HeLa cells using the Abcam ChIP kit protocol. Cells were fixed with formaldehyde for 10 min. ChIP was performed with 2 µg of ab8898 (blue). No antibody was added to the beads control (yellow). Immunoprecipitated DNA was quantified by real time PCR (Taqman approach for active/inactive loci, Sybr green approach for heterochromatic loci). Primers and probes are located in the first kb of transcribed region.



ChIP - ChIP Kit (ab500)

Chromatin immunoprecipitation using ab500 ChIP Kit and Histone H3 (tri methyl K4) antibody (ab12209). Chromatin was prepared from HeLa cells using the Abcam ChIP kit protocol. Cells were fixed with formaldehyde for 10 min. ChIP was performed with 5 µg of ab12209 (blue). No antibody was added to the beads control (yellow). Immunoprecipitated DNA was quantified by real time PCR (Taqman approach for active/inactive loci). Primers and probes are located in the first kb of the transcribed region.



Chromatin immunoprecipitation using ab500 ChIP Kit and Histone H3 antibody (ab1791).

Chromatin was prepared from HeLa cells using the Abcam ChIP kit protocol. Cells were fixed with formaldehyde for 10 min. ChIP was performed with 2 µg of ab1791 (blue). No antibody was added to the beads control (yellow). Immunoprecipitated DNA was quantified by real time PCR (Taqman approach for active/inactive loci, Sybr green approach for heterochromatic loci). Primers and probes are located in the first kb of the transcribed region.

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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