

Product datasheet

ChIP Kit ab500

★★★★☆ 1 Abreviews 35 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
種交差性	交差種: Mouse, Rat, Human
製品の概要	The Protein A ChIP Kit is compatible with rabbit immunoglobulin, various mouse immunoglobulin subtypes, and immunoglobulins from many other species. See the table in the protocol for more details.
アプリケーション	適用あり: ChIP

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

内容	1 kit
Protein A Beads Unblocked	1 x 960µl
Proteinase K	1 x 30µl

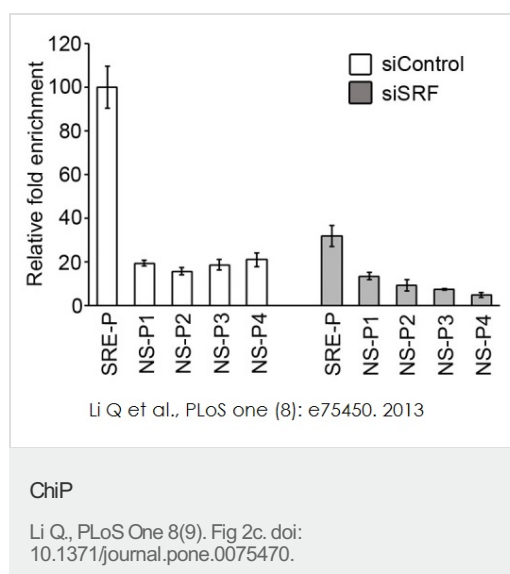
アプリケーション

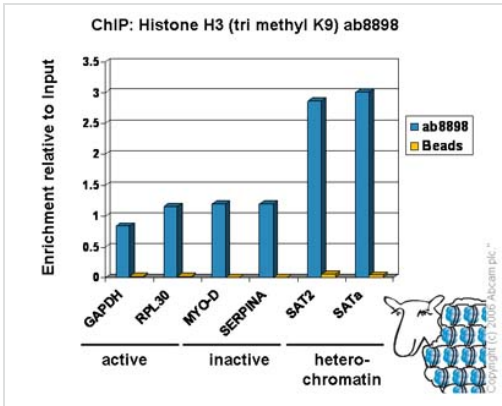
Our [Abpromise guarantee](#) covers the use of **ab500** in the following tested applications.

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.

画像

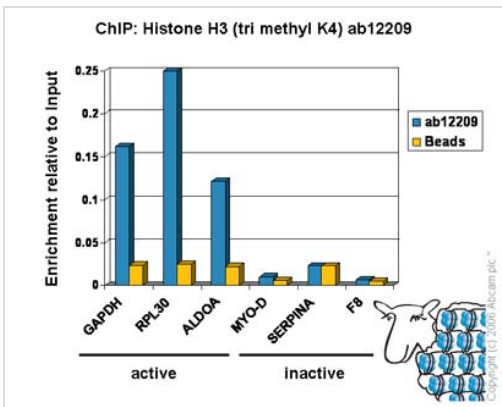




ChIP - ChIP Kit (ab500)

Chromatin immunoprecipitation using ab500 ChIP Kit and Histone H3 (tri methyl K9) antibody (ab8898).

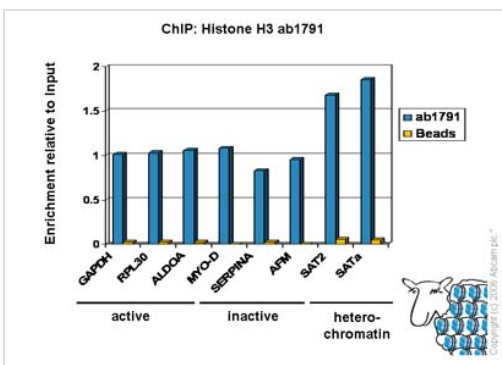
Chromatin was prepared from HeLa cells using the Abcam ChIP kit protocol. Cells were fixed with formaldehyde for 10 min. ChIP was performed with 2 µg of ab8898 (blue). No antibody was added to the beads control (yellow). Immunoprecipitated DNA was quantified by real time PCR (Taqman approach for active/inactive 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.



ChIP - ChIP Kit (ab500)

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.

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 <http://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