

Immunoprecipitation kit ab206996

45 References [画像数 3](#)

製品の概要

製品名	Immunoprecipitation kit
サンプルの種類	Tissue, Adherent cells, Suspension cells
製品の概要	<p>Abcam's Immunoprecipitation Kit (ab206996) can be used to perform immunoprecipitation (IP) and Co-IP for functional studies of immunoprecipitated proteins/complexes and SDS-PAGE or western blot analysis of immunoprecipitated proteins and complexes.</p> <p>Abcam's Immunoprecipitation Kit provides optimized buffers for preparing cell/tissue extracts, antigen binding and washing steps. The Protein A/G Sepharose[®] beads provided in the kit have a higher binding capacity with broader antibody isotype binding than traditional Protein A or Protein G resins.</p>
特記事項	This product is manufactured by BioVision, an Abcam company and was previously called K286 Immunoprecipitation (IP) Kit. K286-25 is the same size as the 25 test size of ab206996.
アプリケーション	適用あり: IP

製品の特性

保存方法 Store at +4°C. Please refer to protocols.

内容	25 tests
10X Wash Buffer X	1 x 20ml
Non-Denaturing Lysis Buffer	1 x 40ml
Protease Inhibitor Cocktail I	1 vial
Protein A/G Sepharose Beads I	1 x 1ml
RIPA Lysis Buffer	1 x 40ml

アプリケーション

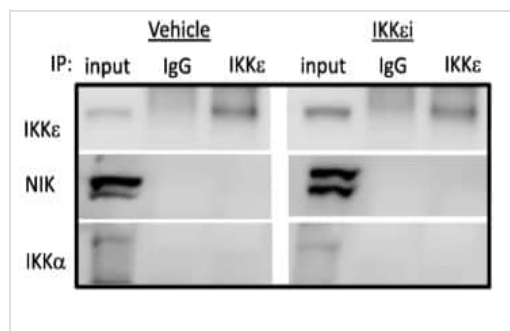
The Abpromise guarantee

Abpromise保証は、次のテスト済みアプリケーションにおけるab206996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
IP		Use at an assay dependent concentration.

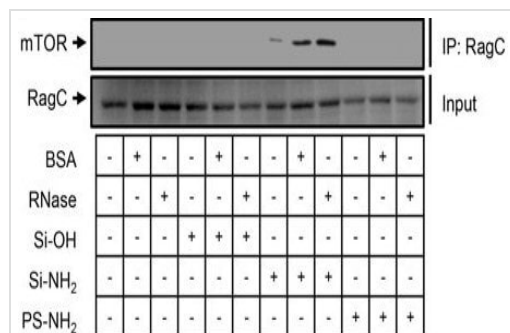
画像



Immunoprecipitation - Immunoprecipitation kit
(ab206996)

Image from House et al., BMC Cancer, 18(1):595; doi: 10.1186/s12885-018-4507-2. Reproduced under the Creative Commons license <https://creativecommons.org/licenses/by/4.0/>

Co-immunoprecipitation experiments indicate IKKε does not interact with NIK or IKKα. Immunoprecipitation was performed on MDA MB 468 cells using the Abcam immunoprecipitation kit (ab206996), according to manufacturer's instructions. Briefly, non-denaturing lysis buffer was used to collect 300 µg of cell lysate was incubated overnight with 3 µg/ml of either control rabbit IgG or IKKε rabbit polyclonal antibody (**ab7891**). Antibody bound proteins were captured using protein A/G sepharose beads, eluted, and analyzed via western blot.

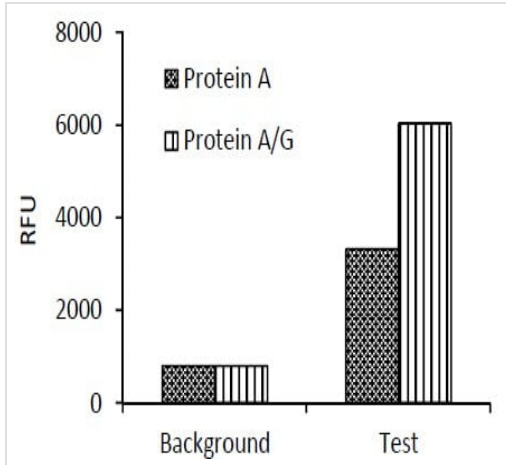


Lunova, Mariia et al.
Scientific reports vol. 7,1 16049. (2017)

Functional Studies - Immunoprecipitation kit
(ab206996)

Lunova, Mariia et al., Scientific reports vol. 7,1 16049., Fig 8, doi:10.1038/s41598-017-16447-6

Huh7 cultured with Si-OH, Si-NH₂, or PS-NH₂ NPs (all 50 µg/ml) bearing BSA or RNase (both 50 µM) as hard protein corona or bare NPs for 4 h. After 4 h post NP treatment, cells were lysed with lysis buffer from immunoprecipitation kit (Abcam). RagC-mTOR complexes were co-immunoprecipitated from the precleared cell lysates with appropriate Ab as described in the manufacturer's instruction. The resulting protein complex was eluted from the beads with Laemmli protein sample buffer for SDS-PAGE and resolved on SDS-PAGE with specific antibody against mTOR



Comparison of Binding Capacity of Protein-A and Protein-A/G Beads.

Comparison of immunoprecipitation using Protein A beads and Immunoprecipitation kit ab206996. HDAC2 activity assay of the antigen-antibody complex captured using Protein A beads or ab206996 by following the same protocol demonstrates that ab206996 is more efficient in IP than Protein A beads.

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