

# Magnetic Conjugation Kit ab269890

2 References 画像数 5

医薬用外毒物

### 製品の概要

製品名	Magnetic Conjugation Kit
製品の概要	<p>Magnetic Conjugation Kit (ab269890) allows conjugation of antibodies to specially treated magnetic particles without the need for extensive optimisation of the conjugation reaction.</p> <p>The 0.5 µm specially treated iron oxide magnetic particles offer you the right balance between high surface area and fast separation. The binding capacity of our magnetic particles is greater than other magnetic particles on the market, and twice that of leading competitors. This allows you to reach the desired sensitivity with fewer particles and therefore minimize the risk of nonspecific binding.</p> <p>Covalent attachment of antibodies to magnetic particles is ideal for immunoprecipitation experiments and provides a more stable alternative to noncovalent attachment of antibodies using protein A or protein G resin.</p> <p><b>Features and Benefits:</b></p> <p>Quick and easy to use – save time, no special knowledge required</p> <p>High binding capacity – achieve high sensitivity with less material</p> <p>High coupling efficiency – No antibody loss during conjugation</p> <p>Covalent bond – Highly stable conjugates</p> <p>No extensive optimisation required – Save precious antibody</p> <p>Scalable – Easy transfer from R&amp;D to manufacturing</p> <p>Freeze dried – Long shelf life</p>

### Specifications

**Core:** Iron oxide

**Mean diameter:** 0.5 µm

**Binding capacity:** ≥70 µg of Goat anti-rabbit antibody per mg of beads

## 特記事項

This product is manufactured by Expedeon, an Abcam company. It was previously called Magnetic Conjugation Kit. Expedeon product code 1300-0005 (3 reaction mini kit) is the same as the 3 x 20 µg size of this product, code 1300-0010 (10 reaction mini kit) is the same as the 10 x 20 µg size, and code 1300-0015 (1 reaction midi kit) is the same as the 1 x 200 µg size of this product.

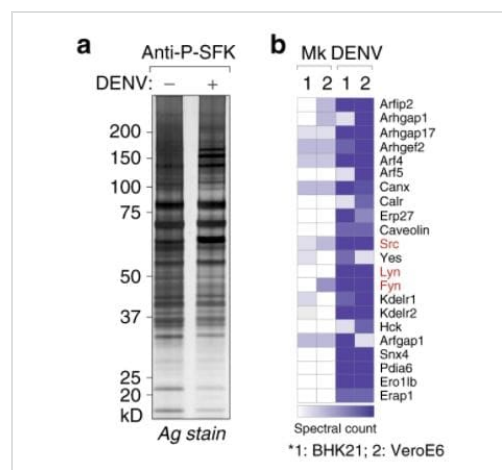
## 製品の特性

### 保存方法

Store at -20°C. Please refer to protocols.

内容	3 x 20 µg	10 x 20 µg	1 x 200 µg
ab273924 - Mag Antibody Diluent	1 x 1500µl	1 x 4ml	1 x 4ml
ab273922 - Mag Quencher	1 x 1500µl	1 x 3ml	1 x 3ml
ab273921 - Mag Storage Buffer	1 x 5ml	1 x 15ml	1 x 15ml
ab273923 - Mag Wash Buffer	1 x 1500µl	1 x 5ml	1 x 5ml
ab273925 - Magnetic particles vial	3 x 20µg	10 x 20µg	1 x 200µg

## 画像

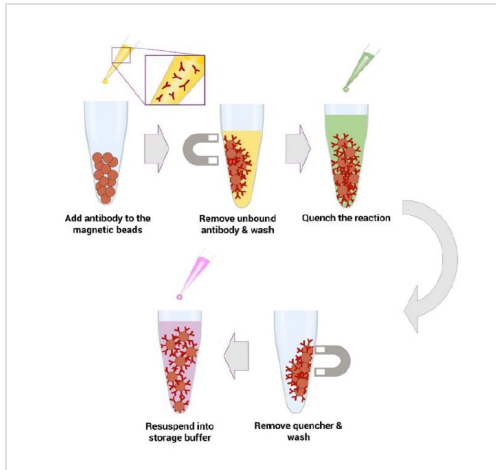


### Mass Spectrometry - Magnetic Conjugation Kit (ab269890)

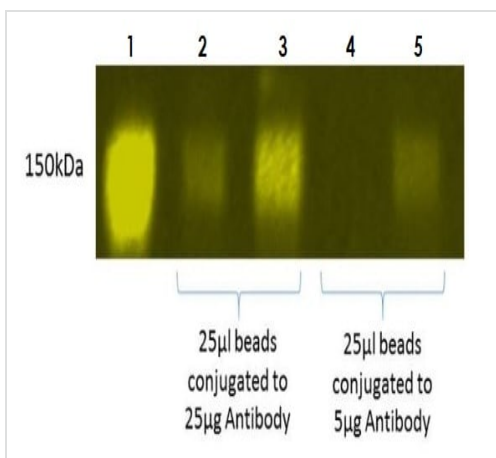
Image from Li et al., Nat Commun., 11(1):5189; doi: 10.1038/s41467-020-19028-w. Reproduced under the Creative Commons license <https://creativecommons.org/licenses/by/4.0/>

Li, Ming Yuan, et al used Magnetic Conjugation Kit (ab269890) as part of screening for SFKs activated upon Dengue infection. They used the kit to conjugate Magnetic beads to anti-phospho-SFK antibodies for use in mass spectrometry.

(a) Large-scale immunoprecipitation of activated SFKs was performed on anti-phospho-SFK antibodies conjugated to the magnetic beads from mock- and Dengue-infected BHK21 and Vero E6 cells (MOI 2, 24 h). Isolated proteins were resolved by SDS-PAGE and detected by silver staining. (b) Entire lanes on gels were sliced into 2-mm sections and subjected to trypsin digest. The peptide mix was processed and analysed by an LTQ Orbitrap mass spectrometer. This allowed the identification of three members Src, Fyn and Lyn, and several SFK-regulated substrates in Dengue-infected samples.

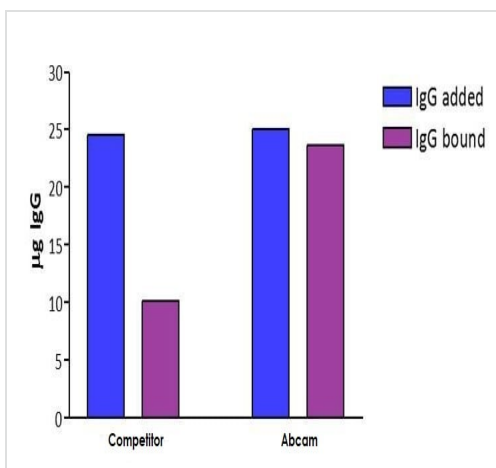


How The Kit Works



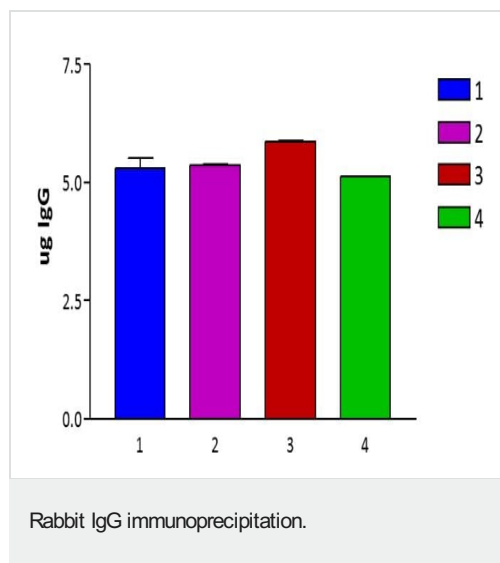
Rabbit IgG immunoprecipitation and detection by Western blot.

25  $\mu$ g and 5  $\mu$ g Goat anti-rabbit antibody were added to each of two Mini vials of magnetic particles, and to 25  $\mu$ L of 1% magnetic beads from a leading competitor, and processed as per the respective protocols. The conjugates were used to immunoprecipitate rabbit IgG from rabbit serum. The eluted samples were run by SDS-PAGE and analysed by Western blot against rabbit IgG. Abcam's Magnetic Conjugation Kit has a greater binding capacity and coupling efficiency than the competitor products, and allows you to detect the antigen of interest with fewer reagents. Lane 1: Rabbit IgG control; Lanes 2 and 4: Competitor; Lanes 3 and 5: Magnetic Conjugation Kit (ab269890).



Magnetic Conjugation Kit binding capacity.

Goat anti-rabbit antibody was conjugated to 25  $\mu$ L of 1% magnetic particles from Magnetic Conjugation Kit (ab269890) (1h procedure), and to magnetic particles from a leading competitor (4h procedure) as per the respective protocols. The quantity of IgG added to the magnetic particle Abcam's Magnetic Conjugation Kit demonstrates a considerably greater binding capacity.



Goat anti-rabbit was labeled with four different batches of magnetic beads. The conjugates were tested in parallel on a Rabbit IgG immunoprecipitation. The bound Rabbit IgG was eluted and quantified by 280nm absorbance.

**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