

# Antibody Concentration And Clean-Up Kit ab102778

**5 References**   [画像数 1](#)

### 製品の概要

---

製品名	Antibody Concentration And Clean-Up Kit
製品の概要	<p>Antibody Concentration And Clean-Up Kit ab102778 allows for the quick and easy concentration of antibodies and proteins using a microcentrifuge-based method. The kit can also be used to reduce the concentration of many unwanted additives often found in antibody formulations such as azide, glycine or Tris.</p> <p>Antibodies are sometimes only available at low concentrations and often contain low molecular weight substances that interfere in labeling reactions with enzymes, biotin, streptavidin and fluorophores. ab102778 can be used to avoid the impact of this.</p> <p>The Antibody Concentration And Clean-Up Kit method utilizes a simple spin column to easily and quickly remove excess buffer from the antibody thereby providing a more concentrated antibody solution. The Antibody Concentration And Clean-Up Kit also allows the experimenter to perform a simple buffer exchange to transfer the antibody into a more favorable buffer for conjugation.</p> <p>Antibodies concentrated using the Antibody Concentration And Clean-up Kit are fully compatible with our Lightning-Link® Antibody Conjugation Kits and our Oligonucleotide Conjugation Kit. However they are not compatible with our nanoparticle conjugation kits (Gold, Latex, Europium and Magnetic).</p> <p>To concentrate antibodies for use with our Gold and Magnetic Conjugation Kits, please use Antibody Concentration And Clean-Up Kit – Gold And Magnetic Nanoparticles (<a href="#">ab204911</a>). To concentrate antibodies for use with our Latex and Europium conjugation kits, please use Antibody Concentration and Clean Up Kit for Latex &amp; Europium (<a href="#">ab269965</a>).</p>

特記事項	This product is manufactured by Expedeon, an Abcam company, and was previously called AbSelect™ Antibody Concentration and Clean-Up Kit. 861-0010 is the same as the 3 units size.
------	--

### 製品の特性

---

## 保存方法

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

内容	3 tests
ab275797 - Conjugation Buffer	1 x 15ml
ab273914 - Spin Cartridge / Collecting Tube Assemblies	3 units

## 画像



Representative Antibody Concentration and Clean-Up Kit (3 columns) (ab102778).

Antibody Concentration And Clean-Up Kit (3 columns) (ab102778)

**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