

## Product datasheet

# APC Conjugation Kit ab201807

1 Image

### Overview

**Product name**

APC Conjugation Kit

**Product overview**

Abcam APC Conjugation Kit (ab201807) uses a simple and quick process to conjugate an antibody to APC. It can also be used to conjugate other proteins or peptides.

To conjugate an antibody to APC using this kit:

- add modifier to antibody and incubate for 3 hours
- add quencher and incubate for 30 mins

The conjugated antibody can be used immediately in WB, ELISA, IHC etc. No further purification is required and 100% of the antibody is recovered for use.

Learn about buffer compatibility below; for incompatible buffers and low antibody concentrations, use our rapid [antibody purification and concentration kits](#). Use the [FAQ](#) to learn more about the technology, or about conjugating other proteins and peptides to APC.

**Notes**

**Amount and volume of antibody for conjugation to APC**

<i>Kit size</i>	<i>Recommended</i>	<i>Maximum antibody</i>
<i>maximum amount of antibody volume</i>		
30 µg	3 x 10 µg	3 x 10 µL
300 µg	3 x 100 µg	3 x 100 µL
1 mg	1 mg	1 mL

**Buffer Requirements for Conjugation**

Buffer should be pH 6.5-8.5.

**Compatible buffer constituents**

If a concentration is shown, then the constituent should be no more than the concentration shown. If several constituents are close to the limit of acceptable concentration, then this can inhibit conjugation.

50mM / 0.6% Tris <sup>1</sup>	0.1%/1% BSA <sup>2</sup>	50% glycerol
0.1% sodium azide	PBS	Potassium phosphate
Sodium chloride	HEPES	Sucrose
Sodium citrate	EDTA	Trehalose

<sup>1</sup> Tris buffered saline is almost always ≤ 50 mM / 0.6%

<sup>2</sup> 1% BSA gives lower quality conjugates, BSA can also interfere with the use of the conjugated antibody in tissue staining.

### ***Incompatible buffer constituents***

ThiomersalProclin Glycine

Arginine GlutathioneDTT

If a constituent of the buffer containing your antibody or protein is not listed above, please check the [FAQ](#) or [contact us](#).

Only purified antibodies are suitable for use, ie. where other proteins, peptides, or amino acids are not present: antibodies in ascites fluid, serum or hybridoma culture media are incompatible.

### **Tested applications**

**Suitable for:** Conjugation

### **Properties**

### **Storage instructions**

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

<b>Components</b>	<b>100 µg</b>	<b>300 µg</b>	<b>30 µg</b>	<b>1 mg</b>
APC Conjugation Mix	1 vial	3 vials	3 vials	1 vial
APC Modifier Reagent	1 vial	1 vial	1 vial	1 vial
APC Quencher Reagent	1 vial	1 vial	1 vial	1 vial

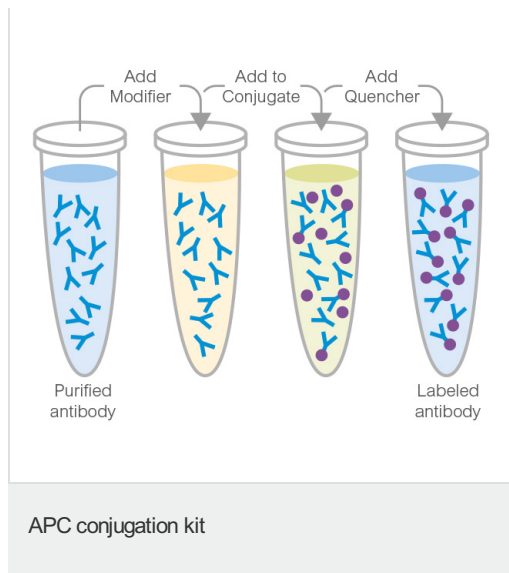
### **Applications**

Our [Abpromise guarantee](#) covers the use of **ab201807** in the following tested applications.

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

<b>Application</b>	<b>Abreviews</b>	<b>Notes</b>
Conjugation		Use at an assay dependent concentration.

### **Images**



**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.com/abpromise> or contact our technical team.

### Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors