

FITC Conjugation Kit ab102884

★★★★★ 2 Abreviews 8 References 画像数 1

製品の概要

製品名

FITC Conjugation Kit

製品の概要

Note: For a faster protocol and the same high quality conjugation, we recommend FITC Conjugation Kit [ab188285](#).

Abcam's FITC Conjugation Kit provides a simple and quick process to conjugate your primary antibodies with FITC. FITC is a green fluorescent dye which has an excitation wavelength of 495nm and an emission wavelength of 519nm.

The conjugated antibody can be used straight away in WB, ELISA, IHC, Immunofluorescence (ICC/IF), Flow cytometry and FRET etc

Learn more about buffer compatibility, protein/secondary antibody conjugation and labelling chemistry in our [FAQs](#).

特記事項

The antibody to be labelled should be purified, in an appropriate buffer for conjugation and at a suitable concentration, as described in the protocol booklet. If not, consider using our [antibody purification and concentration kits](#).

Pre-conjugation considerations

The purified antibody to be labeled should ideally be in 10-50mM amine-free buffer (e.g. MES, MOPS, HEPES, PBS), pH range 6.5 to 8.5. If the buffer is more concentrated or outside this pH range contact the technical support team.

Common non-buffering salts (e.g. sodium chloride), chelating agents (e.g. EDTA), and sugars have no effect on conjugation efficiency. Azide (0.02 to 0.1%) and BSA (0.1 to 0.5%) have little or no effect. Glycerol up to 50% has no effect.

Avoid buffer components that are nucleophilic, as these may react with the kit chemicals. Primary amines (e.g. amino acids or ethanolamine) and thiols (e.g. mercaptoethanol or DTT) fall within this class (Note: Tris-based buffers should be avoided). If your buffer contains primary amines and/or thiols, you should consider using our Concentration and Purification Kits ([ab102778](#) or [ab102784](#)).

Recommended amount and volume of antibody for optimal results

Pack size Amount of antibody Volume of antibody

Up to 20 µg	10-20 µg	4-10 µl
Up to 200 µg	100-200 µg	40-100 µl
Up to 2 mg	1-2 mg	400-1000 µl

Antibody concentrations of 1-4 mg/ml generally give optimal results. For conditions outside the suggested ranges contact the technical support team.

Conjugation reaction

Before adding the antibody to the FITC mix, add 1 µl of Modifier reagent to each 10 µl of antibody to be labelled. Mix gently.

Remove cap from vial of FITC mix and pipette the antibody sample (with added Modifier) directly onto the lyophilized material. Resuspend *gently* by withdrawing and re-dispensing the liquid once or twice using a pipette.

Replace cap on the vial and leave standing in the dark at room temperature (20-25°C) for 3 hours. Conjugations can also be set up and left overnight; longer incubation times have no negative effect on the conjugate.

After incubating for 3 hours (or more), add 1 µl of Quencher reagent for every 10 µl of antibody used and mix gently. The conjugate can be used after 30 minutes. The conjugates do not require purification.

Storage of conjugates

Storage at 4°C is recommended for any conjugate. A preservative may be desirable for long-term storage. Other storage conditions may also be satisfactory. The best conditions for any particular conjugate must be determined by experimentation.

Disclaimer

Labeling of the antibody will not work if the conjugation blocks the active paratope. This situation is rare.

アプリケーション

適用あり: Conjugation

製品の特徴

保存方法

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

内容	300 µg	30 µg	1 mg
FITC mix	3 x 100µg	3 vials	1 vial
Modifier reagent	1 vial	1 vial	1 vial
Quencher reagent	1 vial	1 vial	1 vial

アプリケーション

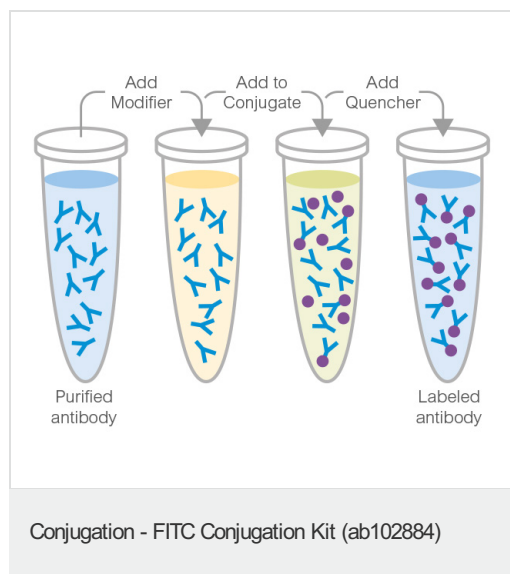
Our [Abpromise guarantee](#) covers the use of **ab102884** in the following tested applications.

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

Conjugation

Use at an assay dependent dilution.

画像



This illustration demonstrates a general procedure and will slightly vary dependent on the conjugate used.

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 <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