

# Streptavidin Atto 542 Monovalent Antibody Labeling Kit ab272188

画像数 3

### 製品の概要

---

**製品名** Streptavidin Atto 542 Monovalent Antibody Labeling Kit

**製品の概要** Streptavidin Atto 542 Monovalent Antibody Labeling Kit ab272188 is used to label biotinylated antibodies with Atto 542.

Atto 542 is very bright and can be used as an alternative to Cy3<sup>®</sup>.

The Antibody - Streptavidin Atto 542 complex produced with this kit is stable and can be used in the same way as antibodies conjugated directly with fluorochromes.

Several antibodies can be labeled with this kit, and its associated kits, and used together in the same experiment.

#### Use this kit to:

- label biotinylated antibodies with fluorescent dyes, and use them in immunostaining just like a directly dye conjugated antibody

- label two or more biotinylated antibodies with different fluorescent dyes\* and then use them together in the same experiment

\*multiple kits required, see list below

- label as little as 0.1 µg of biotinylated antibody in less than 5 mins without washes

#### Technical notes

The 50 µg size of this kit is typically sufficient to label 50 µg of antibody, the 25 µg size is sufficient to label 25 µg of antibody.

When the antibody - streptavidin complex is produced with this kit, due to the monovalent nature of the streptavidin and the blocking solution, there is no risk of forming aggregates of multiple antibodies.

The antibody - streptavidin complex is also very stable, if 2 or more biotinylated antibodies are

labeled with different dyes, and then mixed, no dye will transfer between the antibodies.

### List of available monovalent streptavidin antibody labeling kits

Streptavidin Alexa Fluor® 488 Equivalent ([ab272187](#))

Streptavidin Atto 542 (ab272188) – this kit

Streptavidin Alexa Fluor® 594 ([ab272189](#))

Streptavidin Alexa Fluor® 647 Equivalent ([ab272190](#))

Note: the streptavidin in this kit contains very small amounts of non-monovalent streptavidin, when used with the blocking buffer included in the kit, this does not affect the conjugated antibodies.

#### アプリケーション

**適用あり:** Flow Cyt, IHC

#### 製品の特性

##### 保存方法

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

内容	50 µg	25 µg
Biotin block buffer	1 x 500µl	1 x 250µl
Streptavidin Atto 542 labeling reagent (Monovalent)	1 x 500µl	1 x 250µl

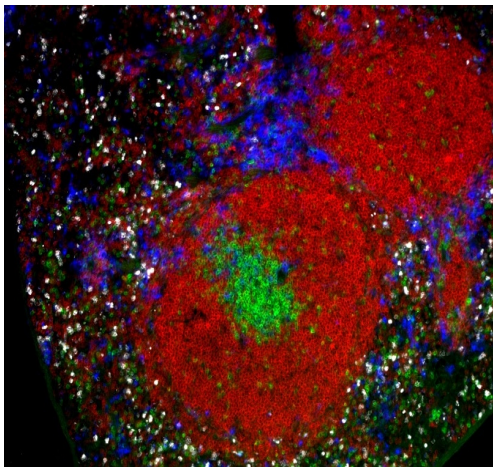
#### アプリケーション

**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ab272188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。

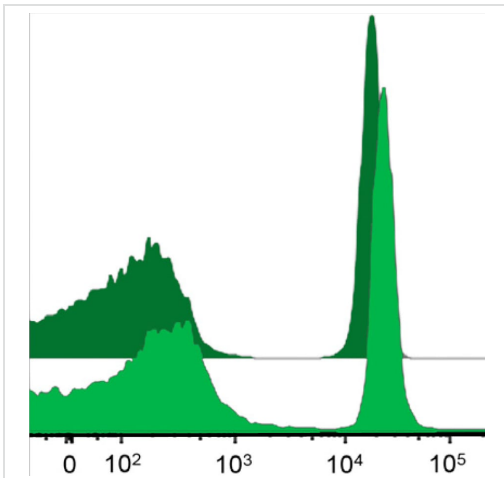
アプリケーション	Abreviews	特記事項
Flow Cyt		Use at an assay dependent concentration.
IHC		Use at an assay dependent concentration.

#### 画像



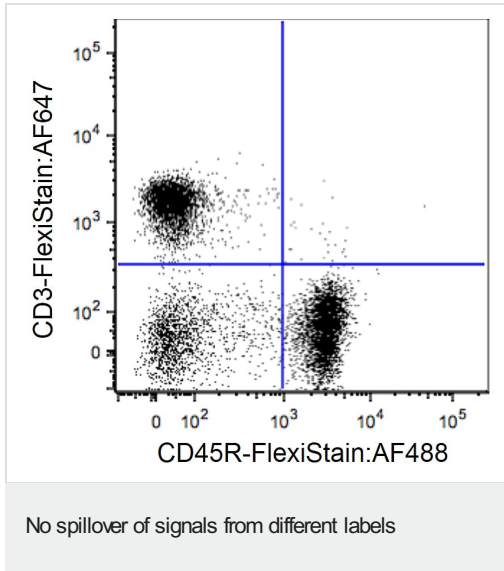
Multiplex 4 color staining

Multiplex immunostaining of a frozen tissue section of mouse spleen using four biotinylated antibodies labeled using Streptavidin (Monovalent) kits. T cells (rat anti-CD3, 7  $\mu\text{g}/\text{ml}$ , green), B cells (rat anti-CD45R, 7  $\mu\text{g}/\text{ml}$ , red), macrophages (rat anti-CD11b, 7  $\mu\text{g}/\text{ml}$ , white), and dendritic cells (hamster anti-CD11c, 7  $\mu\text{g}/\text{ml}$ , blue) were detected using Streptavidin (Monovalent) kits with four different fluorochromes. Picture was taken with a standard fluorescence microscope equipped with appropriate filters for optimal multicolor separation.



Brightness compared to directly conjugated antibodies

Antibodies labeled with Streptavidin (Monovalent) kits are as bright as directly conjugated antibodies. Flow cytometry analysis of splenic B lymphocytes immunolabeled with antibodies against CD45R (B220, clone RA3-6B2, 7  $\mu\text{g}/\text{ml}$ ). Dark green = directly conjugated antibody, light green = Streptavidin (Monovalent)-labeled antibody.



No spillover of signals between Streptavidin (Monovalent)-labeled antibodies. Flow cytometry analysis of T and B cells immunolabeled (7 µg/ml) with Streptavidin Alexa Fluor® 647 (Monovalent) (**ab272190**) and Streptavidin Alexa Fluor® 488 (Monovalent) (**ab272187**).

**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