

Antibody Biotinylation Check Kit ab270030

画像数 2

製品の概要

製品名	Antibody Biotinylation Check Kit
製品の概要	<p>Antibody Biotinylation Check Kit (ab270030) is a quick immunochromatography test that allows you to confirm the successful antibody biotinylation in one easy step, and without the need for any specialized or costly equipment. Unlike traditional methods of assessing the success of biotinylation, such as HABA assays, which can be tedious and time consuming, this kit requires only 10 minutes to perform and very small quantities of your antibody conjugate.</p> <p>The key component of the kit is a nitrocellulose membrane containing a Test line of immobilized Protein A and Protein G with high affinity for the Fc region of a variety of IgG molecules, and a 40 nm Gold-Streptavidin conjugate for detection. When the biotinylated antibody is mixed with the Gold-Streptavidin conjugate, it forms a complex that flows along the nitrocellulose membrane and binds to the Protein A and Protein G on the Test line. A visible red line appears on the strip confirming successful conjugation.</p> <p>Benefits:</p> <p>Quick and easy qualitative test to check success of conjugation</p> <p>Compatible with antibody conjugates generated with Lightning-Link® Biotin conjugation kits (ab201795 and ab201796)</p> <p>No costly equipment required</p> <p>Small volumes of conjugate required (20 µl)</p> <p>General notes:</p> <p>This product is manufactured by Expedeon, an Abcam company. It was previously called Biotin Check&Go! Antibody Labeling Confirmation Kit. Expedeon product code 4001-0030 is the same as the 30-test size of this product.</p>

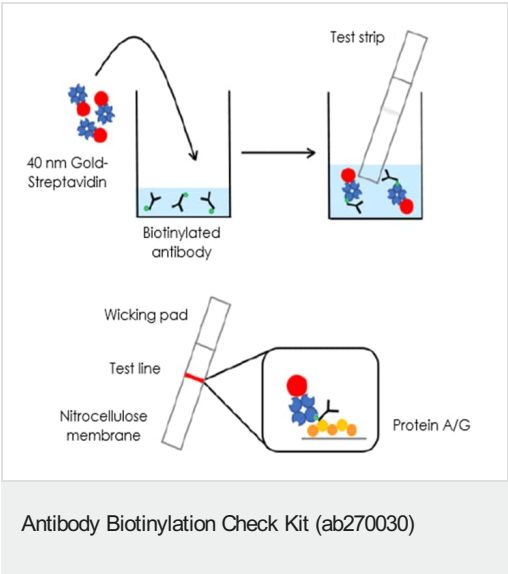
製品の特性

保存方法

Please refer to protocols.

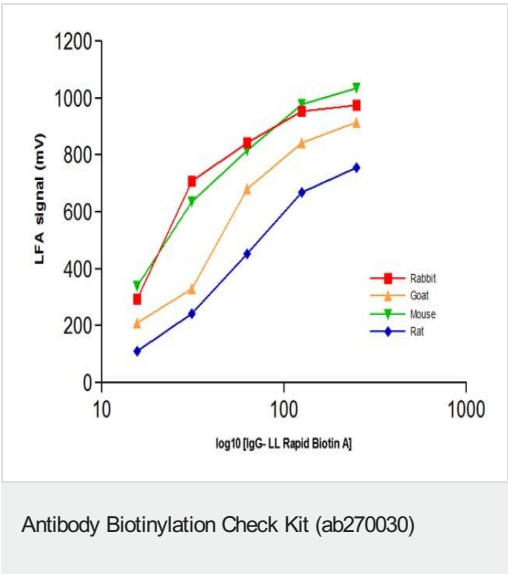
内容	30 tests
ab274084 - 40nm gold conjugated streptavidin (Lyophilized)	1 x 100µg
ab274085 - Biotinylated Rabbit IgG (Lyophilized)	1 x 100µg
ab274081 - Protein A/G Strips	1 x 30 tests
ab274082 - Running Buffer (10X)	2 x 1500µl

画像



How the kit works:

The nitrocellulose membrane contains a Test line of immobilized Protein A and Protein G. Both Protein A and Protein G have a high affinity for the Fc region of a variety of IgG molecules. The membrane also contains an absorbent pad to promote and control the flow of sample through the nitrocellulose. The biotinylated antibody is run on the Protein A/G Strip together with 40 nm gold conjugated streptavidin as the detection label. Biotin binds to streptavidin with exceptionally high affinity and when the biotinylated antibody flows along the nitrocellulose, upon binding to the Protein A and Protein G concentrated on the Test line, a visible red line appears on the strip.



Polyclonal antibodies from rabbit, goat, mouse and rat have been conjugated with Lightning-Link® Rapid Biotin A and run on Protein A / G strips at different concentrations, together with 0.4OD gold conjugated streptavidin. Antibodies from all species provided a red visible line on the strips.

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