

Version 1 Last updated 14 July 2020

ab270051

Streptavidin Antibody Conjugation Check Kit

A product of Expedeon, an
Abcam company

Applicable to Expedeon product codes: 4004-0030

View ab270051

Streptavidin Antibody Conjugation Check Kit:

www.abcam.com/ab270051

(use www.abcam.cn/ab270051 for China, or www.abcam.co.jp/ab270051 for Japan)

For the confirmation of the success of antibody biotinylation in one easy step. Suitable for use with Lightning-Link® Streptavidin antibody labeling kit.

This product is for research use only and is not intended for diagnostic use.

Table of Contents

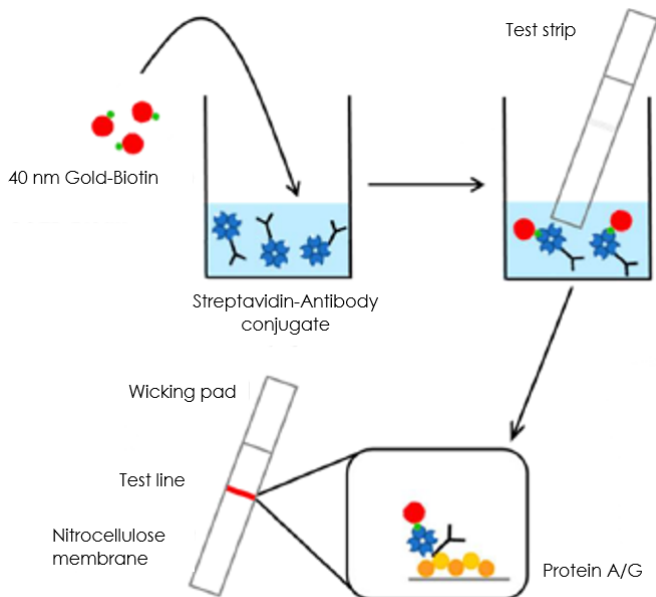
1. Overview	2
2. Materials Supplied and Storage	2
3. Technical Considerations	4
4. Assay Procedure	5

1. Overview

Streptavidin Antibody Conjugation Check Kit (ab270051) is a quick and simple immunochromatography test that allows you to confirm the successful conjugation of streptavidin to your antibody without the need for any specialized or costly equipment.

The key component of the kit is a nitrocellulose membrane containing a 'Test line' of immobilized Protein A and Protein G called a "half strip". Both Protein A and Protein G have a high affinity for the Fc region of a variety of IgG molecules. The "half strips" also contain an absorbent pad to promote and control the flow of sample through the nitrocellulose. The streptavidin-antibody conjugate is run on the Protein A/G strip together with 40 nm gold conjugated Biotin as detection label. Biotin binds streptavidin with exceptionally high affinity and when the streptavidin-antibody conjugate flows along the nitrocellulose membrane, upon binding to the Protein A and Protein G on the T-line, a visible red line appears on the strip.

Streptavidin Antibody Conjugation Check Kit (ab270051) is suitable for use with Lightning-Link® Streptavidin antibody labelling kit.



2. Materials Supplied and Storage

Store Protein A/G Strips at +4°C. Store all other components at -20°C immediately on receipt. Kit can be stored for 1 year from receipt, if components have not been reconstituted.

Avoid repeated freeze-thaws of reagents.

Item	30 tests	Storage temperature
10X Running Buffer	2 vials	-20°C
Positive Control – Streptavidin-Rabbit IgG Conjugate (lyophilized)	1 vial	-20°C
40 nm Gold Biotin (lyophilized)	1 vial	-20°C
Protein A/G Strips	30 Strips	+4°C

Not supplied: 96-wells low binding plate, Bovine serum albumin (BSA)

3. Technical Considerations

Protein A and Protein G affinity for immunoglobulins:

Species	Ig subclasses	Binding to Protein A	Binding to Protein G
Rabbit	IgG	High	High
Human	IgG ₁	High	High
	IgG ₂	High	High
	IgG ₃	No affinity	High
	IgG ₄	High	High
	IgA	Low	No affinity
	IgD	Low	No affinity
	IgE	Low	No affinity
Human	IgM	Low	No affinity
Pig	IgG	High	High
Mouse	IgG ₁	Low/Medium	Medium
	IgG _{2a}	High	High
	IgG _{2b}	High	Medium
	IgG ₃	Low/Medium	Medium
	IgM	Low	No affinity
Goat	IgG	Low	High
Sheep	IgG	Low	High
Rat	IgG	Low	High
	IgG ₁	Low	Low/Medium
	IgG _{2a}	Low	High
	IgG _{2b}	Low	Low/Medium
	IgG _{2c}	Low	Low/Medium
	IgM	Low	No affinity

Δ Note: Low/no affinity for a specific IgG subclass may lead to low/no signal but the conjugate may be fine.

4. Assay Procedure

- Equilibrate all materials and prepared reagents to room temperature just prior to use and gently agitate.
- 4.1** Dilute the 10x Running Buffer with distilled water down to 1x
- 4.2** Add 0.1% BSA (final concentration) to 1x Running Buffer
- 4.3** Reconstitute the 40 nm Gold Biotin with 150 μL of 1x Running Buffer + 0.1% BSA to obtain 2 OD
- 4.4** Dilute your sample(s) in 1x Running Buffer + 0.1% BSA; you will need 10 μL for each test; the optimal conjugate concentration range is 250 - 15.6 ng/mL. It is recommended to run the samples in duplicate and include the positive control (see below).
- 4.5** For a single test mix 10 μL of diluted streptavidin-antibody, 4 μL of reconstituted Gold Biotin 2 OD and 26 μL of 1x Running Buffer + 0.1% BSA and mix through pipetting.
- 4.6** Load 40 μL /well of your mix in a 96-well non-sticky plate or a suitable container.
- 4.7** Dip the strip in the well.
- 4.8** Run for 10 minutes. Check the result by eye. A red line will become visible if the conjugation was successful.

Positive control – Streptavidin-Rabbit IgG Conjugate

- 4.9 Reconstitute the vial in 300µl of 1x Running Buffer + 0.1% BSA
- 4.10 For a single run mix 10 µL of positive control, 4µL of reconstituted Gold Biotin 2 OD and 26 µL of 1x Running Buffer + 0.1% BSA and mix through pipetting.
- 4.11 Load 40 µL/well of positive control.
- 4.12 Dip the strip in the well.
- 4.13 Run for 10 minutes.

The positive control will produce a visible red line indicating a successful dipstick assay.

Technical Support

Copyright © 2020 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

Austria

wissenschaftlicherdienst@abcam.com | 019-288-259

France

supportscientifique@abcam.com | 01.46.94.62.96

Germany

wissenschaftlicherdienst@abcam.com | 030-896-779-154

Spain

soportecientifico@abcam.com | 91-114-65-60

Switzerland

technical@abcam.com

Deutsch: 043-501-64-24 | Français: 061-500-05-30

UK, EU and ROW

technical@abcam.com | +44(0)1223-696000

Canada

ca.technical@abcam.com | 877-749-8807

US and Latin America

us.technical@abcam.com | 888-772-2226

Asia Pacific

hk.technical@abcam.com | (852) 2603-6823

China

cn.technical@abcam.com | 400 921 0189 | +86 21 2070 0500

Japan

technical@abcam.co.jp | +81-(0)3-6231-0940

Singapore

sg.technical@abcam.com | 800 188-5244

Australia

au.technical@abcam.com | +61-(0)3-8652-1450

New Zealand

nz.technical@abc.com | +64-(0)9-909-7829