

ab270029 Streptavidin Gold Conjugate (20 nm, 100D)

A product of Expedeon, an
Abcam company

Applicable to Expedeon product codes 251-0200, 251-1000.

View ab270029

Streptavidin Gold Conjugate (20 nm, 100D) datasheet:

www.abcam.com/ab270029

(use www.abcam.cn/ab270029 for China, or www.abcam.co.jp/ab270029 for Japan)

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	3
3. Formulation	3
4. Instructions	3

1. Overview

Abcam's Streptavidin Gold Conjugate (20 nm, 100D) is a unique coated nanoparticle that has an avidly bound protective surface coat which can withstand the most extreme conditions (e.g. 2.5M NaOH at 70°C for >1 hour).

By covalently attaching streptavidin to the surface coat the streptavidin ligand is irreversibly attached and cannot desorb, which is a major problem with products in which the streptavidin is passively adsorbed to the metal.

Streptavidin Gold Conjugate binds with high affinity to biotinylated molecules.

2. Materials Supplied and Storage

Store product at +4°C upon receipt.

Item	Quantity		Storage temperature
Streptavidin Gold Conjugate	200 µL	1 mL	+4°C

3. Formulation

Gold Streptavidin is shipped in a buffer composed of TBS containing 0.1% detergent. If you wish to exchange the Gold Streptavidin into a specific buffer for your assay or test, centrifuge the conjugate at 9,000 x *g* during 20 minutes. Carefully remove the supernatant and add your preferred buffer. It is important to avoid substances that have a very high affinity for gold (e.g. thiols).

The maximum absorbances for the 20 nm Gold Streptavidin is at 528 nm. To determine the effective concentration of the conjugate obtained we advise to measure the Absmax of light using an UV-vis spectrophotometer after diluting your sample to an appropriate range for your piece of equipment.

4. Instructions

A common application for 20 nm Streptavidin Gold Conjugates is to be used as a detector in lateral flow assays (LFAs). They can also be exploited as a LFA control, when biotinylated BSA (available separately) is immobilized on the C-line of a strip.

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.

For all technical or commercial enquiries please go to:

www.abcam.com/contactus

www.abcam.cn/contactus (China)

www.abcam.co.jp/contactus (Japan)