



# InnovaCoat® GOLD Streptavidin 10nm, 20nm and 40nm 10 OD

Applicable to:

250-0200	40nm 200µl pack
251-0200	20nm 200µl pack
252-0200	10nm 200µl pack

250-1000	40nm 1ml pack
251-1000	20nm 1ml pack
252-1000	10nm 1ml pack

Release 7

20/03/2017

## Introduction

At the core of this product is InnovaCoat® GOLD, 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.

InnovaCoat® GOLD Streptavidin binds with high affinity to biotinylated molecules.

## Shipping and storage conditions

InnovaCoat® GOLD Streptavidin is shipped at ambient temperature in a tamper-evident polypropylene container.

On receipt, store the product at +4°C. The shelf life is >12 months.

## Formulation

InnovaCoat® GOLD Streptavidin is shipped in a buffer composed of TBS containing 0.1% detergent.

If you wish to exchange the InnovaCoat® GOLD Streptavidin into a specific buffer for your assay or test, centrifuge the conjugate according to the table below:

Gold nanoparticles size	Centrifuge speed	Time
10nm	21,500g	45 minutes
20nm	9,000g	20 minutes
40nm	9,000g	6 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 10nm, 20nm and 40nm IC-GOLD Streptavidin are at 520nm, 528nm and 530nm respectively. To determine the effective concentration of the conjugate obtained we advise to measure the Abs<sub>max</sub> of light using an UV-vis spectrophotometer after diluting your sample to an appropriate range for your piece of equipment (e.g. if a 40nm gold conjugate is at 10 OD and is diluted 1:10 the Abs<sub>530nm</sub> for a 1 cm light path is expected to be around 1 AU).

## Limited use license

Innova Biosciences' products are offered for research purposes alone, and are not intended for human, therapeutic or diagnostic use. The purchase of this product conveys to the buyer (whether the buyer is a not-for-profit, academic or for-profit entity) the non-transferable right to use the amount of product purchased and the components of the product for in-house research. The buyer shall not sell or otherwise transfer this product, its components, or materials prepared therefrom to any third party. The buyer shall not use this product or its components for commercial purposes. For the avoidance of doubt, 'commercial purposes' means any activity by a party for consideration and includes, without limitation, use of the product or its components (i) in the manufacturing of other materials (ii) to provide a service, information or data, (iii) for therapeutic, diagnostic or prophylactic purposes, or (iv) for repackaging/resale, whether or not such product or its components are resold for use in research. The use of this product by the buyer constitutes agreement with the terms of this limited use label.

Many of Innova Biosciences' products and technologies are protected by patents or are the subject of patent filings. For information on purchasing a license for commercial use please contact our Business Development Office, Babraham Hall, Babraham, Cambridge, UK, CB22 3AT. Tel +44(0)1223 661000; e-mail info@innovabiosciences.com

## Instructions

A common application for 20nm and 40nm 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.

The 10nm Streptavidin gold conjugates are more suitable for imaging in electron microscopy.

For additional applications please contact us at: <https://www.innovabiosciences.com/contact-us.html>

## Related products and services

- InnovaCoat® GOLD Biotin 10 OD 40nm (240-0200; 240-1000).
- InnovaCoat® GOLD Custom Services, including optimization of covalent antibody conjugates, covalent attachment of small molecules and ligands.
- Biotinylated BSA (2050-0020)

Visit our website: [www.innovabiosciences.com](http://www.innovabiosciences.com) for more information.