

InnovaCoat® GOLD Goat Anti-Mouse IgG 80nm, 40nm, 20nm and 10nm 10 OD

Applicable to:	217-0200	200µl	217-1000	1ml
	216-0200	200µl	216-1000	1ml
	215-0200	200µl	215-1000	1ml
	214-0200	200µl	214-1000	1ml

Release 1. © EXPEDEON. 22/02/2018

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).

The Goat Anti-Mouse IgG antibody (GAM) is irreversibly attached to the surface coat via a covalent bond. This irreversible attachment doesn't allow the desorption of the antibody, which is a major problem with products in which the antibody is only passively adsorbed to the metal.

The InnovaCoat® GOLD conjugated to GAM antibody recognizes Mouse IgG which is often used as a primary antibody in various research applications. For this reason secondary antibody anti-mouse IgGs can be useful tools for the analysis of target proteins.

SHIPPING AND STORAGE CONDITIONS

InnovaCoat® GOLD GAM 80, 40, 20 and 10nm at 10 OD concentration are shipped at ambient temperature in a small tamper-evident polypropylene container. On receipt, store the product at +4°C. The shelf life is >12 months.

FORMULATION

InnovaCoat® GOLD GAM 80nm, 40nm, 20nm and 10nm 10 OD are shipped in a buffer composed of TBS containing 0.1% detergent. If you wish to exchange the InnovaCoat® GOLD GAM 80, 40, 20 and 10 nm into a specific buffer for your assay or test, centrifuge the conjugate in a microfuge at:

- 9,000g 6 minutes (80 and 40nm GAR GOLD)
- 9,000g 20 minutes (20 nm GAR GOLD)
- 21,500g 45 minutes (10nm GAR GOLD)

Carefully remove the supernatant, gently tap the pellet and add your preferred buffer. It is important to avoid substances that have a very high affinity for gold (e.g. thiols).

INSTRUCTIONS

The InnovaCoat® Gold labelled with GAM is suitable for the following applications:

When using InnovaCoat® Gold GAM we recommend to try several dilutions according to the specific application. If an incubation of InnovaCoat GOLD GAM with the target antibody is performed, do not centrifuge the gold complex after this stage (an antibody titration is recommended).

Product Code	Size	LFA	Flow Cytometry	DLS	SEM	TEM
217-0200	80					
217-1000	nm	•	•	•		
216-0200	40					
216-1000	nm	•	•		•	•
215-0200	20					
215-1000	nm	•	•		•	•
214-0200	10					
214-1000	nm		•		•	•

The table above is provided merely as a guideline. For additional applications please contact our Technical Support.

RELATED PRODUCTS AND SERVICES

Description	Prod. Code
InnovaCoat® GOLD Goat Anti-Rabbit IgG 80nm 10 OD	(221-0200; 221-1000)
InnovaCoat® GOLD Goat Anti-Rabbit IgG 40nm 10 OD	(213-0200; 213-1000)
InnovaCoat® GOLD Goat Anti-Rabbit IgG 20nm 10 OD	(219-0200; 219-1000)
InnovaCoat® GOLD Goat Anti-Rabbit IgG 10nm 10 OD	(218-0200; 218-1000)
InnovaCoat® GOLD Custom Services including optimization of covalent antibody conjugates, covalent attachment of small molecules and ligands.	

TECHNICAL SUPPORT

For technical enquiries get in touch with our technical support team at: technical.enquiries@expedeeon.com
For further information see our website: www.expedeeon.com

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, 25 Norman Way, Over Cambridgeshire CB24 5QE, United Kingdom Tel +44 (0)1223 873364; e-mail info@expedeeon.com