

Colloidal gold nanoparticles 10, 20, 40, 80nm

Applicable to: 3001-0020, 3001-0100, 3010-0020, 3010-0100
 3201-0010, 3201-0100, 3210-0010, 3210-0100, 3215-0010, 3215-0100
 201-0010, 201-0100, 210-0010, 210-0100, 212-0010, 212-0100, 220-0010, 220-0100
 3801-0010, 3801-0100, 3810-0010, 3810-0100, 3815-0010, 3815-0100

Release 2. © EXPEDÉON. 27/11/2018

INTRODUCTION

Expédeon's colloidal gold nanoparticles are ultra-high quality, with a narrow size distribution and uniform shape, supplied in a range of sizes and concentrations for maximum flexibility.

SPECIFICATIONS

Particle size	Effective diameter (TEM)	Standard deviation	% CV
10nm	10nm ± 1nm	< 1nm	< 10%
20nm	20nm ± 2nm	< 2nm	< 10%
40nm	40nm ± 4nm	< 4nm	< 10%
80nm	80nm ± 8nm	< 8nm	< 10%

CONTENTS

Bottles of gold nanoparticles:

10nm

(3001; 3010)-0020 20ml bottle at 1 or 10 OD

(3001; 3010)-0100 100ml bottle at 1 or 10 OD

20nm

(3201; 3210; 3215)-0010 10ml bottle at 1, 10 or 15 OD

(3201; 3210; 3215)-0100 100ml bottle at 1, 10 or 15 OD

40nm

(201; 210; 212; 220)-0010 10ml bottle at 1, 10, 15 or 20 OD

(201; 210; 212; 220)-0100 100ml bottle at 1, 10, 15 or 20 OD

80nm

(3801; 3810; 3815)-0010 10ml bottle at 1, 10 or 15 OD

(3801; 3810; 3815)-0100 100ml bottle at 1, 10 or 15 OD

STORAGE

The colloidal gold nanoparticles are shipped at ambient temperature, upon receipt store between 2-8°C, do not freeze.

RELATED PRODUCTS

InnovaCoat® GOLD nanoparticle conjugation kits provide a simple, easy to use approach to conjugation of antibodies to gold nanoparticles. Details of the full range of InnovaCoat® products can be found on our website.

Description	Code
InnovaCoat® GOLD 10nm nanoparticle conjugation kit	228-
InnovaCoat® GOLD 20nm nanoparticle conjugation kit	229-
InnovaCoat® GOLD 40nm nanoparticle conjugation kit	230-
InnovaCoat® GOLD 80nm nanoparticle conjugation kit	231-

InnovaCoat® GOLD Secondary Antibody conjugates provide a simple and reliable way of targeting primary antibodies for immunological techniques. Details of the full range of InnovaCoat® GOLD Secondary Antibody conjugates products can be found on our website.

InnovaCoat® GOLD Custom Services, including optimization of covalent antibody conjugates, covalent attachment of small molecules and ligands.

Limited use license

Use and Use Restrictions: The Products are sold, and deliverables of any services are provided for the purposes of the buyer's internal in vitro research, development or educational use only, not for in vivo, or any therapeutic or diagnostic use, nor for resale, or for providing services or any other commercial use of any kind, including without limitation, for any transfer in any form (including as part of a kit) to a third party; for analysis or reverse engineering of the Product or for manufacturing. Products should only be used in accordance with any safety data sheets, guidance or protocols that we issue from time to time and are available for download from our Site. Protective clothing should be used at all times when handling our Products. Safety datasheets relating to all Products are available for download from the Site or upon request.

Expédeon grants no other license or rights under any intellectual property in respect of Products or services deliverables and in particular grants no license to use any Product or deliverables for any commercial purposes. Sale of Products or service deliverables by us or our authorised distributors are expressly conditional upon the customer's agreement with these restrictions, which the customer gives upon placing an order for Products or deliverables. If you wish to use any Product or deliverables for any purpose other than your own internal research as described above, you will require an additional licence from Expédeon. Please contact licensing@expedeon.com.