

Gold Conjugation Kit (10 nm, 20 OD) ab201808

[8 References](#) [画像数 2](#)

製品の概要

製品名	Gold Conjugation Kit (10 nm, 20 OD)
製品の概要	<p>Abcam's Gold Conjugation Kit allows antibodies or proteins to be covalently attached to ultra-stable Gold nanoparticles at very high OD quickly and easily. The nanoparticles in the Abcam's Gold Conjugation Kit have a protective surface coat that can withstand the most extreme conditions (e.g. 2.5M NaOH at 70°C for > 1 hour). The hands-on time for the Abcam Gold conjugation procedure is around 2 minutes and the conjugate is ready to use within 20 minutes. The researcher simply pipettes the biomolecule into a vial containing the nanoparticles of the Gold Conjugation Kit.</p> <p>The nanoparticles in this kit are supplied as a freeze-dried mixture. The conjugation reaction is initiated simply by reconstituting the freeze-dried nanoparticles with the antibody, which becomes attached (through lysine residues) to the surface of the nanoparticles.</p> <p>The resulting covalent conjugates are more stable than those prepared by passive adsorption methods. Moreover, unlike passive methods, the coating process is independent of the isoelectric point of the antibody, avoiding the need for extensive trials at different pH values. All antibodies can be labelled at a single pH.</p> <p>Learn more about buffer compatibility, protein/secondary antibody conjugation and labeling chemistry in our FAQs.</p> <p>Benefits:</p> <p>Easy and rapid conjugation – only 2 minutes hands-on time and 100% recovery of materials</p> <p>Site-specific labelling – Antigen binding sites of antibodies are free to bind the target molecule</p> <p>Proprietary surface coating prevents metal-protein interactions, and enables covalent attachment to the Gold – Stable conjugates formed</p> <p>Fully scalable – Easy transfer from R&D to manufacturing</p> <p>Uniform spherical shape and narrow size distribution – Consistent high quality and excellent</p>

batch-to-batch reproducibility

Buffer requirements:

The biomolecule to be conjugated should ideally be in 10 mM amine-free buffer (e.g. MES, MOPS, HEPES), pH range 6.5 to 8.5. Sugars have no effect on conjugation efficiency. For incompatible buffers and low antibody concentrations, use our rapid **antibody purification and concentration kits** for Nanoparticles. To learn more about incompatible buffers, please refer to the protocol booklet.

特記事項

This product is manufactured by Expedeon, an Abcam company, and was previously called 10nm InnovaCoat[®] GOLD. 228-0015 is the same as the 1 x 25 µg size. 228-0010 is the same as the 10 x 2.5 µg size. 228-0005 is the same as the 3 x 2.5 µg size.

The 3 and 10 Test Conjugation Kits are designed to label 12 µl of antibody per vial. The 1 Test Conjugation Kit is designed to label 120 µl of antibody per vial.

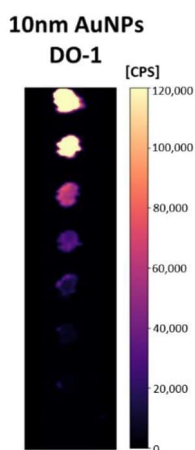
製品の特性

保存方法

Store at -20°C. Please refer to protocols.

内容	3 x 2.5 µg	10 x 2.5 µg	1 x 25 µg
ab273944 - Gold 10nm	3 x 2500ng	10 x 2500ng	1 x 25µg
ab273943 - Gold Antibody Diluent	1 x 1ml	1 x 4ml	1 x 4ml
Gold Quencher Reagent	1 x 700µl	1 x 700µl	1 x 700µl
Gold Reaction Buffer	1 x 750µl	1 x 750µl	1 x 750µl

画像

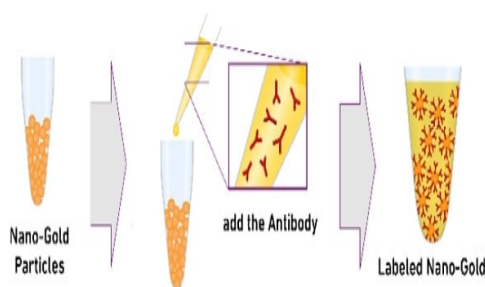


Dot Blot - Gold Conjugation Kit (10 nm, 20 OD)

(ab201808)

Image from Vlcovska, Marcela, et al., *Molecules*, 26(3):630; doi: 10.3390/molecules26030630. Reproduced under the Creative Commons license <https://creativecommons.org/licenses/by/4.0/>

Vlcovska, Marcela, et al used Gold Conjugation Kit (10 nm, 20 OD) (ab201808) as part of examining antibody activity after binding to nanoparticles. They used the kit to conjugate Gold to monoclonal anti-DO-1 antibody for use in LA-ICP-MS dot-blot analysis. Inductively-coupled laser ablation plasma mass spectrometry (LA-ICP-MS) dot-blot analysis of p53 protein standard (total protein amount 0.10-10.00 ng) using DO-1 antibody conjugates with Europium nanoparticles (EuNPs).



GOLD Conjugation Kit (10nm, 20OD)

GOLD Conjugation Kit (10nm, 20OD)

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.co.jp/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors