

Product datasheet

GOLD Conjugation Kit (80nm, 20 OD) ab154876

[1 Image](#)

Overview

Product name GOLD Conjugation Kit (80nm, 20 OD)

Product overview 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 hands-on time for the Gold Conjugation procedure is about 2 minutes and the conjugate is ready to use within 20 minutes.

The Gold nanoparticles in this kit are supplied as a freeze dried mixture. The conjugation reaction is initiated simply by adding a solution of the antibody, which becomes attached (via lysine residues) to the Gold surface.

The resulting covalent conjugates are more stable than those prepared by passive adsorption methods. Moreover, unlike passive methods, the coating procedure is not dependent on the isoelectric point of the antibody, and extensive trials at different pH values are not required; all antibodies react at a single fixed pH.

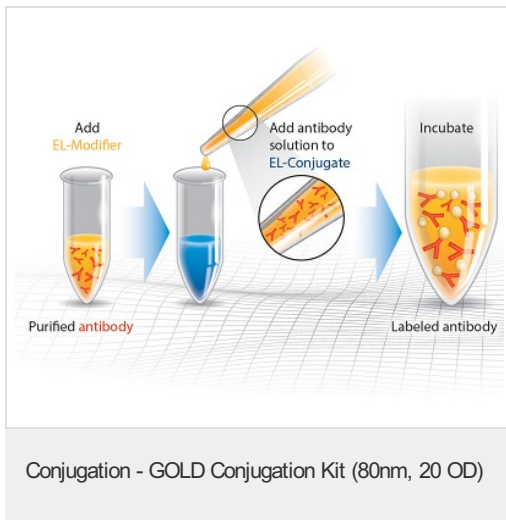
Notes The 3 and 10 Test Conjugation Kits are designed to label 12 µl per vial.
The 1 Test Conjugation Kit is designed to label 120 µl per vial.

Properties

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

Components	3 tests	10 tests	1 tests
Gold	3 vials	10 vials	1 vial
Gold Antibody Diluent	1 vial	1 vial	1 vial
Gold Quencher Reagent	1 vial	1 vial	1 vial
Gold Reaction Buffer	1 vial	1 vial	1 vial

Images



Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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.com/abpromise> or contact our technical team.

Terms and conditions

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