

LANCE® *Ultra*

Eu-W1024 labeled Anti-Human IgG Antibody

Product number: AD0212 **Lot Number:** 3249105

Product Format: AD0212: 10 µg

AD0074: 50 µg

AD0075: 1 mg

Manufacturing date: August 2, 2023 **Document version:** 1

Product Information

Storage Buffer: 50 mM Tris-HCl pH 7.4, 150 mM NaCl, 0.05% NaN₃, 0.1% BSA

Molecular Weight: 160 000

Storage: Store at 4°C.

Stability: This product is stable for at least 30 months from the manufacturing date when stored in its original packaging and the recommended storage conditions.

Quality Control

The QC release specifications are based on spectrophotometric analysis of the labeled antibody. We certify that results meet our quality release criteria.

Labeling Ratio: 6.52 (Eu/protein)
Concentration: 100 µg/mL (0.625 µM)

Please visit our website for additional information on LANCE *Ultra* technology at www.revvity.com

The information provided in this document is valid for the specified lot number and date of analysis. This information is for reference purposes only and does not constitute a warranty or guarantee of the product's suitability for any specific use. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume any liability for any errors or damages arising from the use of this document or the product described herein. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.

AD0074-5-0212-R Rev01