

Research Use Only. Not for use in diagnostic procedures.

Eu-N1 labeled Anti-goat IgG Antibody

Product No.: AD0303, AD0304, AD0305

Material Provided

Format: AD0303, 50 µg
AD0304, 200 µg
AD0305, 1 mg

Manufacturing Date:

Product Information

Antibody: The antibody is a mouse monoclonal anti-goat antibody. It reacts with goat IgGs. Cross-reactivity with human, mouse, rabbit, rat, and bovine immunoglobulins is less than 0.2% and cross reactivity with sheep immunoglobulin is less than 0.5% when determined with DELFLIA.

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

Molecular Weight: 160 000

Stability: This product is stable for at least **24 months** from the manufacturing date when stored in its original packaging and the recommended storage conditions.

Storage Conditions: Store at 4°C

Quality Control

The QC release specifications are based on spectrophotometric analysis and a fluorescence assay performed on a EnVision™ instrument. We certify that these results meet our quality release criteria.

Labeling Ratio: 19-30 (Eu/Ab)

Concentration: 1000 µg/mL (6.25 µM)

This product is not for resale or distribution except by authorized distributors.

LIMITED WARRANTY: PerkinElmer warrants that, at the time of shipment, the above-named product is free from defects in material and workmanship and conforms to the specifications set forth above. PerkinElmer makes no other warranty, express or implied with respect to the product and expressly disclaims any warranty of merchantability or fitness for any particular purpose. Notification of any breach of the foregoing warranty must be made within 60 days of receipt of the product, unless otherwise provided in writing by PerkinElmer. No claim shall be honored if the customer fails to notify PerkinElmer within the period specified. The sole and exclusive remedy of the customer for any breach of the foregoing warranty is limited to either the replacement of the non-conforming product or the refund of the invoice price of the product. PERKINELMER SHALL NOT BE LIABLE FOR ANY DIRECT, INDIRECT, SPECIAL, INCIDENTAL, CONSEQUENTIAL OR PUNITIVE DAMAGES, WHETHER BASED ON CONTRACT, TORT, STRICT LIABILITY OR OTHERWISE, ARISING OUT OF THE DESIGN, MANUFACTURE, SALE, DELIVERY, OR USE OF THE PRODUCTS, EVEN IF THE LIMITED REMEDIES PROVIDED HEREIN FAIL OF THEIR ESSENTIAL PURPOSE OR PERKINELMER IS ADVISED OF THE POSSIBILITY OF SUCH DAMAGES.

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
P: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs