

Research Use Only. Not for use in diagnostic procedures.

DELFLIA Eu-N1 Goat Anti-HRP Antibody

Product No.: AD0306, AD0307

Lot No.: XXX-XXX-X

Material Provided

Format: AD0306 25 µg
AD0307 100 µg

Manufacturing Date: XXXXXXX

Product Information

Antibody: DELFLIA Eu-N1 goat anti-HRP antibody binds strongly to HRP-labeled mouse, rabbit, goat, and human IgGs as well as HRP-labeled streptavidin. It has < 0.25% cross-reactivity with un-labeled human, mouse, rat, rabbit, donkey, goat, sheep, and bovine IgG.

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

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 an EnVision™ instrument. We certify that these results meet our quality release criteria.

Labeling Ratio: XXX (Eu/Ab)

Concentration: 100 µg/mL (0.625 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

Copyright© 2012, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners