

DELFI[®]**Eu-N1-labeled Anti-Phosphotyrosine (PY20) Antibody****Product number:** AD0039 **Lot Number:** 3385251**Product Format:**
AD0038: 50 µg
AD0039: 1.0 mg**Manufacturing date:** 3/24/2025 **Document version:** 1**Product Information**

Antibody: The PY20 antibody is a mouse monoclonal antibody that binds to phosphorylated tyrosine residues.

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

Molecular Weight: 160 000

Storage: Store at -20°C. Repeated freezing and thawing should be avoided.

Stability: This product is stable for at least 28 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 and a fluorescence assay performed on a VICTOR[™] instrument. We certify that these results meet our quality release criteria.

Labeling Ratio: 7.7 (Eu/Ab)
Concentration: 1000 µg/mL (6.25 µM)

Please visit our website for additional information on DELFIA 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.

AD0038-9-R Rev01