

DELFI[®]

Eu-N1-labeled Anti-c-myc Antibody

Product number: **AD0112** Lot Number: 3363084

Product Format: AD0112: 50 µg
AD0113: 1.0 mg

Manufacturing date: **12/19/2024** Document version: 1

Product Information

Antibody: The antibody is a mouse monoclonal antibody. It recognizes the epitope sequence (EQKLISEEDL), which was derived from the human c-myc protein.

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: 8.5 (Eu/Ab)
Concentration: 500µg/mL (3.125 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.

AD0112-3-R Rev01