

Delfia®

EuTDA Cytotoxicity Reagents

Product number: AD0116 Lot Number: 3311298

Product Format: 960 assay points

Expiration date: 06/26/2025 Document version: 1

Product Information

Reagent	Item Number	Quantity
BATDA Reagent	C136-100	50 µL
Lysis Buffer	AD0116-A	0.5 mL
Europium Solution	AD0116-B	200 mL
Microtitration Strips, uncoated	AD0116-C	10 plates

Intended Use: This product is intended for loading of cells to be used in short term cytotoxicity tests performed utilizing time-resolved fluorometry in the detection.

Storage: Store at 4°C

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.

AD0116-R Rev 01