

Delfia®

# EuTDA Cytotoxicity Reagents

Product number: AD0116 Lot Number: 3261469

Product Format: 960 assay points

Expiration date: 10/11/2024 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](http://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