

## BATDA Reagent

**Product Number:** C136-100

**Lot Number:** 3311853

**Production Date:** 17 OCT 2024

---

### Material Provided

**Format:** 50 µL

### Product Information

**Description:** The DELFLIA® BATDA Reagent is intended for labeling of cells to be used in short term cytotoxicity tests performed utilizing time-resolved fluorometry in the detection.

**Storage Buffer:** Dimethylsulphoxide (DMSO)

**Stability:** This product is stable for at least **12 months** from the manufacturing date when stored in its original packaging and the recommended storage conditions.

**Storage Conditions:** Store at 4°C.

### Quality Control

**hydrolysis test** tda-h<sub>2</sub>o : >7  
batda-h<sub>2</sub>o

**cytotoxicity test** max signal : >3.7  
background

The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. 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

C136-100-R REV01