

DELFLIA®

Lysis Buffer

Product number: 4005-0010 Lot Number: 3228353

Product Format: 4005-001: 30 mL

Manufacturing date: 12/06/2023 Document version: 1

Product Information

Description: 1 bottle (30 mL) of Lysis Buffer containing digitonin (0.03 %) and 19 %dimethylsulphoxide (DMSO).

Application: This product is intended for cell lysis procedures in time-resolved fluorometry tests.

Storage: Store at +2 - +8°C.

Stability: This product is stable for at least 21 months from the manufacturing date when stored in its original packaging and the recommended storage conditions.

Reconstitution: Heat up the Lysis Buffer on water bath (+37°C) prior to use. Digitonin in the Lysis Buffer may precipitate during storage. In case of a white precipitate, it is preferable to warm the Lysis Buffer to +50 - +60°C to completely dissolve digitonin. Should precipitation occur, make sure that it is dissolved before use. Let the reagent reach room temperature (+20 - +25°C) before use.

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.

4005-0010-R Rev 01