

AlphaLISA®

# CETSA® Cell Lysis Buffer Selection Kit

**Product number:** CETSA-BUF-KIT **Lot Number:** 3206818

**Kit Contents:** CETSA® Cell Lysis Buffer 1 (5X), 10 mL  
 CETSA® Cell Lysis Buffer 2 (5X), 10 mL  
 CETSA® Cell Lysis Buffer 3 (5X), 10 mL  
 CETSA® Cell Lysis Buffer 4 (5X), 10 mL  
 CETSA® Cell Lysis Buffer 5 (5X), 10 mL

**Manufacturing date:** September 21, 2023 **Document version:** 1

## Product Information

**Application:** The CETSA® Cell Lysis Buffer Selection Kit contains CETSA® Cell Lysis Buffers 1-5 and has been validated for use with Alpha CETSA® and Alpha SureFire® CETSA® assays.

**Storage:** Store the Alpha CETSA® Lysis Buffer Selection Kit, and all of its components at 4 °C.

**Stability:** The undiluted stock solutions of each buffer are stable for at least 12 months from the manufacturing date when stored in original packing at 4 °C.

Please note that CETSA® is a registered trademark of Pelago Bioscience AB who hold the exclusive rights to the CETSA® patent family. If you do not have yet a valid license, please contact Pelago Biosciences at [sales@pelagobio.com](mailto:sales@pelagobio.com) to discuss your needs.

Please visit our website for additional information on Alpha CETSA® 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.

CETSA-BUF-KIT-R Rev01