

## Assay buffer without detergents (5x conc.)

**Product Number:** CR85-100

**Lot Number:** 3281382

**Production Date:** 22 APR 2024

---

This product is provided for DELFLIA® cell based assays or in assays where the presence of detergents may interfere. DELFLIA® Assay Buffer without detergents is five times concentrated.

### PRODUCT CONTENTS:

250mL. It contains 250 mmol/L Tris-HCl, pH 7.8, 4.5% NaCl, 0.25% NaN<sub>3</sub>, 0.25% Bovine-g-globulin, 100µmol/L DTPA, 2.5% BSA and an inert red dye.

### STORAGE

Store @ 2-8 °C. Material should be used within 15 months of the production date.

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

CR85-100-R REV01