

## Stabilizer (DTPA-purified BSA 7.5%)

**Product Number:** CR84-100

**Lot Number:** 3203992

**Production Date:** 06 NOV 2023

---

Gives low background as a blocking reagent in the DELFLIA® assays.

The stabilizer is intended for use in coating of plates and as a stabilizing agent for storage of lanthanide-labeled compounds. Gives low background as a blocking reagent in the DELFLIA® assays. It consists of bovine serum albumin (BSA) purified from trace amounts of heavy metals.

### PRODUCT CONTENTS:

10 nmol in 0.075 ml 50 mM Tris-HCl (pH 7.8) 0.9% NaCl with <0.1% sodium azide as a preservative and 0.1% BSA

### 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

CR84-100-R REV01