

## DELFLIA<sup>®</sup> Inducer 4013-0010

### INTENDED USE

The DELFLIA<sup>®</sup> Inducer is an acidic chelating detergent solution intended for use in the rapid quantitative determination of Eu<sup>3+</sup>/Sm<sup>3+</sup> in time-resolved fluoroimmunoassays.

The DELFLIA Inducer is suitable for use in multilabel assays that require DELFLIA Enhancer (prod. no. C500-100) for development of Tb<sup>3+</sup>/ Dy<sup>3+</sup> fluorescence.

The solution dissociates Eu<sup>3+</sup>/Sm<sup>3+</sup> from solid phase bound Eu/Sm-labeled antibodies during a time period of a few minutes to form a homogeneous and highly fluorescent Eu/Sm-(BFPP)<sub>3</sub>(TOPO)<sub>2-3</sub> micellar chelate solution. The solution allows highly sensitive Eu<sup>3+</sup>/Sm<sup>3+</sup> measurements to be performed when using a time-resolved fluorometer.

### PACKAGE CONTENTS

1 bottle (250 mL) of DELFLIA Inducer

The DELFLIA Inducer is a ready-for-use solution with Triton X-100<sup>1</sup>, glycine, hydrochloric acid and chelators.

Store at +2 - +8°C until expiry date stated on the bottle label. Shelf life 2 weeks in dark, at room temperature (+20 - +25°C). **Protect from light.**

### PRODUCT USE

Allow the bottle to reach room temperature before use.

Avoid contamination of any kind by minimizing contact with the solution.

Use e.g. an "Eppendorf Multipette" type of pipette (prod. no. 1296-014) with 5 mL Combitips (prod. no. 1296-016), or alternatively the DELFLIA Plate Dispense (prod. no. 1296-041). If other dispensing systems are used, tubing that comes in contact with the DELFLIA Inducer should be made of teflon.

After incubation with Eu-labeled antibodies, wash the solid phase with neutral washing solution (saline). Add 200 µL of DELFLIA Inducer directly from the reagent bottle to each well. **When automatic dispensers are used make sure that the tubing is thoroughly flushed with fresh DELFLIA Inducer before dispensing into the well.**

The DELFLIA Inducer is intended to be used with the N1, DTPA, W1024 and W2014 lanthanide chelates.

---

DELFLIA is a registered trademark of PerkinElmer, Inc.

<sup>1</sup> Triton is a registered trademark of the Rohm and Haas Co.

When using these DELFIA lanthanide chelate labeled reagents, shake the frame slowly for 5 minutes on the DELFIA Plateshake (prod. no. 1296-003) using slow shaking (950 rpm). Then measure the fluorescence with a time-resolved fluorometer. The fluorescence is stable for several hours if evaporation is prevented. However, we recommend measurement within 1 hour as external factors may cause a decrease in signal with time, although this is extremely rare.

In case of DELFIA Inducer evaporation, the fluorescence signal can be returned by dispensing DELFIA Inducer into the well, but the signal level may be higher than after first dispensing.

The same Combitip must not be used for pipetting any other reagent.

When using the 1420 VICTOR™ Multilabel Counter<sup>2</sup> which has not been normalized previously with DELFIA Inducer-based reagents, run Eu/Sm Dual Label Normalization by using DELFIA Inducer-based reagents before dual label measurement (Eu<sup>3+</sup> and Sm<sup>3+</sup>).

**NOTE:** Normalization will overwrite the existing normalization data. Prepare a new plate for DELFIA Inducer-based reagents. Make a copy of existing plate and run Eu/Sm normalization with DELFIA Inducer-based reagents.

## PRODUCT SPECIFICATIONS

The maximum Eu<sup>3+</sup> background in this solution is  $3 \times 10^{-13}$  mol/L.

April 2002



---

<sup>2</sup> VICTOR is a trademark of PerkinElmer, Inc.