

## DELFLIA<sup>®</sup> Enhancement Solution 4001-0010

The DELFLIA<sup>®</sup> Enhancement Solution is an acidic chelating detergent solution intended for use in the quantitative determination of  $\text{Eu}^{3+}/\text{Sm}^{3+}$  when using time-resolved fluoroimmunoassay. The solution dissociates  $\text{Eu}^{3+}/\text{Sm}^{3+}$  from solid phase bound Eu-labelled antibodies during a time period of a few minutes to form a homogeneous and highly fluorescent  $\text{Eu}/\text{Sm}-(2\text{-NTA})_3(\text{TOPO})_{2-3}$  micellar chelate solution. The solution allows highly sensitive  $\text{Eu}^{3+}/\text{Sm}^{3+}$  measurements to be made when using the time-resolved fluorometer. The DELFLIA Enhancement Solution is also used as part of  $\text{Tb}^{3+}/\text{Dy}^{3+}$  measurement together with DELFLIA Enhancer (prod. no. C500-100).

### PACKAGE CONTENTS

1 bottle (1000 mL) of Enhancement Solution

The Enhancement Solution is a ready for use solution with Triton X-100<sup>1</sup>, acetic acid and chelators.

Store at +2 - +8°C until expiry date stated on the bottle label. Shelf life 6 months at room temperature (+20 - +25°C). Protect from light when not in regular use.

### 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 Multipipette" 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, Teflon tubings are preferred in cases where the Enhancement Solution comes in contact with the tubing.

After incubation with Eu-labelled antibodies, wash the solid phase with neutral washing solution (saline). Add 200  $\mu\text{L}$  of Enhancement Solution directly from the reagent bottle to each well. When automatic dispensers are used make sure that the tubing is thoroughly flushed with fresh Enhancement Solution before dispensing into the well.

Shake the frame slowly for 5 minutes on the DELFLIA Plateshake (prod. no. 1296-003) using slow shaking (950 rpm). Then measure the fluorescence with the 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.

DELFLIA is a registered trademark of Wallac Oy.

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



The Enhancement Solution should be dispensed using only the recommended dispensers. The same Combitip must not be used for pipetting any other reagent.

## SYMBOLS



For *in vitro* diagnostic use



Expiry date



Store at



Lot no.



Manufacturer