

DELFIA[®] Samarium Standard Solution B115-100

INTENDED USE

The Samarium Standard Solution is intended for *in vitro* use in the adjustment of Wallac time-resolved fluorometers.

PACKAGE CONTENTS

1 bottle (50 mL) of 10 nmol/L Samarium Standard Solution. The concentration of 10 nmol/L Samarium Standard solution is 10.0 ± 0.6 nmol/L.

The Samarium Standard Solution is a solution where samarium is in a fluorescent form in the Enhancement Solution.

Store **in dark** at +2–+8 °C until expiry date stated on the bottle label. The expiry date is valid only when the product is stored tightly closed and protected from light at +2–+8 °C. Shelf life 3 months at room temperature (+19–+25 °C). To avoid the risk of contamination, use the contents of the bottle immediately after opening.

WARNINGS AND PRECAUTIONS

Disposal of all waste should be in accordance with local regulations.

PRODUCT USE

Allow the bottle to reach room temperature (+19–+25 °C) before use.

Take a clean plate or similar and dispense Samarium Standard Solution to each well with a clean dispenser (not a glass pipette), as specified in the service manual for the instrument. Measure the fluorescence within 30 minutes with the samarium measurement parameter.

The avoidance of samarium contamination and resulting high fluorescent background demands high standard pipetting technique. The same pipette tip must not be used for pipetting any other reagent. In case of contamination, first use Enhancement Solution (prod.no. 1244-105) for cleaning, then wipe with a soft cloth or tissue paper soaked in clean water.

The target signal for adjustment is typically around 100 000 counts, when measured in DELFIA Microtitration Plates (prod. no. C551-110) using the factory preset samarium label parameters.

Last revision December 2009

Manufactured by:

Wallac Oy
Mustionkatu 6
FI-20750 Turku
FINLAND