

DELFLIA[®] Terbium Standard Solution C558-100

INTENDED USE

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

PACKAGE CONTENTS

1 bottle (50 mL) of 1 nmol/L Terbium Standard Solution

The Terbium Standard Solution is a ready for use solution with terbium, cetyltrimethylammonium bromide, Triton X-100¹, acetic acid Tris base, sodium acetate and chelators.

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 at +2 - +8°C. Shelf life 3 months at room temperature (+20 - +25°C).

PRODUCT USE

Allow the bottle to reach room temperature (+20 - +25°C) before use.

Take a clean plate and dispense 250 µL of Terbium Standard Solution to each well with a clean dispenser. Measure the fluorescence within 30 minutes with the terbium measurement parameter.

The avoidance of terbium 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 terbium contamination, firstly use Enhancement Solution (prod. no. 1244-105) for cleaning, then wipe with a soft cloth or tissue paper soaked in clean water.

PRODUCT SPECIFICATION

The signal obtained is typically 500 000 counts, when measured in DELFIA Microtitration Plates (prod. no. C551-110) using the factory preset terbium label parameters.

Last revision March 2005

Manufactured by:

PerkinElmer Life and Analytical Sciences,
Wallac Oy
P.O. Box 10
FIN-20101 Turku
FINLAND