

DELFLIA[®] 1 nmol/L Europium Standard Solution B119-100

INTENDED USE

Europium 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 Europium Standard Solution. The concentration of 1 nmol/L Europium Standard Solution is 1.00 ± 0.08 nmol/L.

Europium Standard Solution is a ready-for-use solution with europium, Triton¹ X-100, acetic acid 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 tightly closed and protected from light at +2 - +8 °C. The shelf life of the product is 3 months at room temperature (+19 - +25 °C). To avoid the risk of contamination, use the contents of the bottle immediately after opening.

INTRODUCTION

Europium Standard Solution is a solution where europium is in a fluorescent form in the Enhancement Solution.

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 Europium 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 factory preset europium measurement parameter.

The target signal for the adjustment is 1 000 000 (\pm 80 000) cps when measured in DELFLIA Microtitration Plates (prod. no. C551-110) using the factory preset europium label parameters.

The avoidance of europium 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 for cleaning, then wipe with a soft cloth or tissue paper soaked in clean water.

Last revision May 2009

Manufactured by:

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