

DELFLIA[®] Assay Buffer 4002-0010

For Research Use Only

INTENDED USE

The DELFLIA[®] Assay Buffer is a buffered protein and detergent solution intended for use in dissociation-enhanced time-resolved fluoroimmunoassays (DELFLIA) that include Eu/Sm/Tb-labeled antibodies or antigens. It is optimized to give a minimum non-specific background in solid phase assays.

PACKAGE CONTENTS

1 bottle (1000 mL) of Assay Buffer

Assay Buffer is ready for use Tris-HCl buffered NaCl solution (pH 7.8), containing < 0.1 % NaN₃, bovine serum albumin (BSA), bovine gamma globulins, Tween 40, diethylenetriaminepentaacetic acid (DTPA), and an inert red dye.

Store at +2 - +8°C until expiry date stated on the bottle label.

PRODUCT USE

The DELFLIA Assay Buffer is meant for use as a diluent for Eu/Sm/Tb-labeled compounds. Labeled reagents are diluted to the required final concentration with the Assay Buffer just prior to the immunoreaction procedure.

Do not store labeled reagents in the Assay Buffer.

SYMBOLS

Expiry date



Store at



Lot no.



Manufacturer

Last revision September 2005

Manufactured by:

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