

DELFLIA[®]

Assay Buffer 1244-106 / 1244-111

INTENDED USE

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

Assay Buffer is available in two volumes:

- 1 bottle of 50 mL (prod. no. 1244-106)
- 1 bottle of 250 mL (prod. no. 1244-111)

Assay Buffer is a 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

For Research Use Only.

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.

Last revision November 2007

Manufactured by:

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

Tel. int. + 358-2-2678 111

DELFLIA is a registered trademark of PerkinElmer, Inc.

