

DELFLIA[®] Streptavidin Microtitration Strips 4009-0010

For Research Use Only

HIGH CAPACITY

INTENDED USE

The DELFLIA[®] Streptavidin Microtitration Strips are intended for use as a solid phase in dissociation-enhanced time-resolved fluorometric assays, e.g. fluoroimmunoassays.

PACKAGE CONTENTS

10 plates (8 x 12 wells) of Streptavidin Microtitration Strips

The wells are directly coated with streptavidin allowing high binding capacity.

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

PRODUCT USE

1. Transfer the required number of microtitration strips to a strip frame. Do not take more strips than can be easily handled within 30 minutes.

Note: Open the foil from three sides only and fold it aside leaving the plate-specific information on the package. Return the remaining strips into the package and press the foil cover back on as tightly as possible. Leave the desiccant in the package. Alternatively, store the remaining strips in a resealable plastic bag with the desiccant.

2. Proceed according to your own optimized assay protocol.

WARNINGS AND PRECAUTIONS

The DELFLIA Streptavidin Microtitration Strips are intended for research use only.

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

WARRANTY

Purchase of this reagent gives the purchaser the right to use this material in his own research. Further distribution of this reagent is expressly prohibited. Purchase of this product implies agreement with these conditions of sale.

Last revision April 2005

DELFLIA is a registered trademark of PerkinElmer, Inc.

Manufactured by:

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

