

# **DELFLIA<sup>®</sup> Anti-rabbit IgG Microtitration Strips 4008-0010**

## **For Research Use Only**

### **INTENDED USE**

The DELFLIA<sup>®</sup> Anti-rabbit IgG Microtitration Strips are intended for use as a solid phase in dissociation-enhanced time-resolved fluoroimmunoassays.

### **PACKAGE CONTENTS**

10 plates (8 x 12 wells) of Anti-rabbit IgG Microtitration Strips

The wells are coated with goat anti-rabbit IgG antibodies.

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 Anti-rabbit IgG 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

Manufactured by:

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