

DELFLIA®

## Eu-labeled Streptavidin

Product number: 1244-360 Lot Number: 3265247

Eu/SA Ratio: 7.4

Expiration date: 03 APR 2025 Document version: 2

### Product Information

#### Intended Use:

DELFLIA® Eu-Labeled Streptavidin is intended for use as a generic reagent to detect biotin-labeled compounds, e.g. antibodies or antigens in various types of immunoassay (1,2) or nucleic acid probes in hybridization assays (3,4) based on solid-phase separation. The label, Eu<sup>3+</sup>, is measured using dissociation-enhanced time-resolved fluorometry (5,6).

#### Product Description:

Streptavidin is derived from *Streptomyces avidinii* and is labeled with the Eu-chelate of N<sup>1</sup>-(p-isothiocyanatobenzyl)-diethylenetriamine-N<sup>1</sup>,N<sup>2</sup>,N<sup>3</sup>,N<sup>3</sup>-tetraacetic acid.

Binding capacity is about 10 µg biotin per mg streptavidin (7).

In contrast to avidin, streptavidin has an isoelectric point close to neutrality and does not contain carbohydrates. Therefore, non-specific binding is less than with avidin.

#### Stability/Storage:

Store at +2 - +8°C.

Please visit our website for additional information on DELFLIA technology at [www.revvity.com](http://www.revvity.com)

The information provided in this document is valid for the specified lot number and date of analysis. This information is for reference purposes only and does not constitute a warranty or guarantee of the product's suitability for any specific use. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume any liability for any errors or damages arising from the use of this document or the product described herein. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.

1244-360-R Rev02