

Lance®

Eu-W8044 labeled streptavidin

Product number: AD0061 **Lot Number:** 3228847

Product Format: 1.0 mg

Expiration date: 07/16/2025 **Document version:** 1

Product Information

Labeling Yield: 6.5
Storage Buffer: 50 mM Tris-HCl buffered saline (pH 7.8) with azide as an antimicrobial agent and 0.1% BSA
Concentration: 500 ug/mL
Labeling Reagent: Lance Terpy Eu DTA chelate
Labeling Conditions: +4°C, pH 9, o/n
Purification Method: Gel filtration
Storage: 2-8 °C (36-46 °F)

Please visit our website for additional information on Lance technology at 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.

AD0061-R Rev01