

AlphaScreen®

Streptavidin Donor Beads

Product number: 6760002B **Lot Number:** 3267378

Material provided: Donor Beads at 5 mg/mL in in 25 mM Hepes, 100 mM NaCl, 0.05% Kathon, pH 7.4

Product Format: 6760002S: 1 mg*, packaged as 1 x 200 µL

6760002: 5 mg, packaged as 1 x 1 mL

6760002B: 50 mg, packaged as 2 x 5 mL

* Note: 1mg translates into 2,000 assay points based on a final bead concentration of 20 µg/mL in a 25 µL/well reaction volume.

Manufacturing date: February 22, 2024 **Document version:** 1

Product Information

Stability: This product is stable for at least 24 months from the Manufacturing date if used and stored under recommended conditions.

Storage: Store undiluted at 4°C protected from light. Freeze-thaw is not recommended and can cause the beads to form aggregates.

Recommended Use: AlphaScreen® Donor beads are light-sensitive. All Alpha assays using the Donor beads should be performed under subdued laboratory lighting (< 100 lux). Green filters (LEE 090 filters can be applied to light fixtures).

Please visit our website for additional information on AlphaLISA 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.

6760002X-R Rev01