

AlphaLISA®

## Acceptor Beads

**Product number:** 6772001, 6772002, 6772003 **Lot Number:** 3190851

**Material provided:** AlphaLISA Acceptor Beads at 20 mg/mL in 100 mM Tris, 0.05% Kathon, pH 8.0

**Product Format:** 6772001: 1 mg\*, packaged as 1 x 50 µL  
6772002: 5 mg, packaged as 1 x 250 µL  
6772003: 50 mg, packaged as 5 x 500 µL

**Manufacturing date:** 10/02/2023 **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.

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

### Quality Control

Europium content of beads is determined and normalized by dividing the Fluorescence of beads at 612 nm by their Absorbance at 525 nm. This ratio is known to be proportional to expected signal intensity of conjugated beads\*. We certify that these results meet our requirements.

*\*Note: Assay performance, including signal intensity, is dependent on the assay configuration. For better performance, we recommend that the Alpha assay be optimized for each assay configuration.*

**TEST:** FU615/A525

**RESULTS:** 1.4

**SPECIFICATION:** ≥ 1.0

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.

677200X-R Rev01