

AlphaPlex™

545 Tb Acceptor Beads

Product number: 6782001 **Lot Number:** 3335707

Material provided: Unconjugated 545 Acceptor Beads at 20 mg/mL in 0.1M MES, pH 6.0

Product Format: 6782001: 1 mg*, packaged as 1 x 50 µL
 6782002: 5 mg, packaged as 1 x 250 µL
 6782003: 50 mg, packaged as 5 x 500 µL

Manufacturing date: September 10, 2024 **Document version:** 1

Product Information

Stability: This product is stable for at least 12 months from manufacturing date if used and stored under recommended conditions.

Storage: Store undiluted at 4°C protected from light. Do not freeze.

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.

678200X-R Rev01