

AlphaPlex™-545

Streptavidin Acceptor Beads

Product number: AP125TB-M **Lot Number:** 3280668

Material provided: AlphaPlex-545 Streptavidin Acceptor Beads at 5 mg/mL in PBS pH 7.2 supplemented with 0.05% Kathon as a preservative.

Product Format: AP125TB-C: 250 µg, 50 µL, 500 assay points

AP125TB-M: 5 mg, 1 mL, 10 000 assay points

AP125TB-R: 25 mg, 5 mL, 50 000 assay points

The number of assay points is based on an assay volume of 25 µL in 384-well assay plates using a final bead concentration of 20 µg/mL.

Manufacturing date: March 27, 2024 **Document version:** 1

Product Information

Application: This product is designed for use as a tool to generate Alpha assays involving biotinylated molecules bound to an acceptor bead.

Storage: Store product in the dark at 4 °C.

Stability: This kit is stable for at least 4 months from the date of manufacture when stored in its original packaging and the recommended storage conditions.

Quality Control

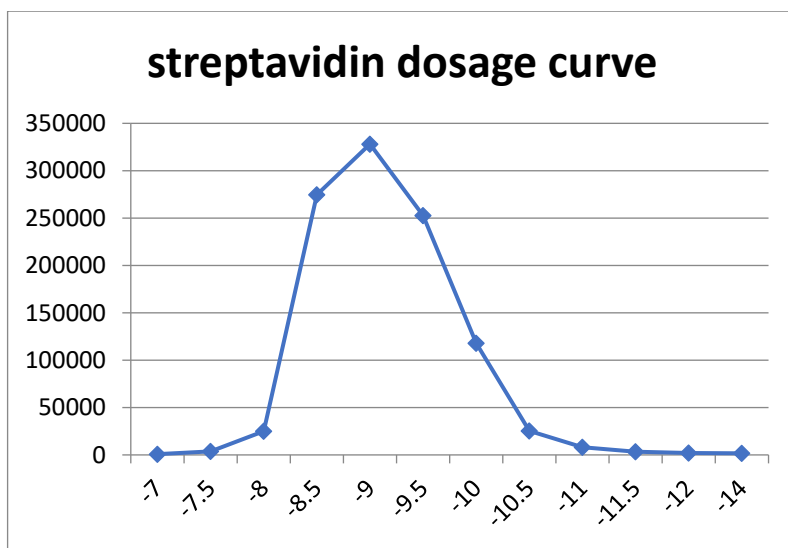
Lot to lot consistency is confirmed in an Alpha assay. Maximum and minimum signals and EC50 were measured on the EnVision Multilabel Plate Reader with Alpha option. We certify that these results meet our quality release criteria. Maximum counts may vary between bead lots and the instrument used, with no impact on assay quality

EC₅₀: 0.50 nM
Min counts: 1562 counts
Max counts: 443795 counts

Recommendations

- Sodium azide should not be added to stock solutions or assay components. Final concentrations of sodium azide higher than 0.001 % will decrease the AlphaPlex signal.
- Total signal varies with temperature and incubation time. For consistent results, identical incubation times and temperature should be used for all plates.
- The AlphaPlex signal is detected with an EnVision Multilabel Reader equipped with the ALPHA option using the following settings: Total Measurement Time: 550 ms, Laser 680 nm Excitation Time: 180 ms, Mirror: D640as, Emission Filter: Wavelength 535nm, bandwidth: 40nm, Transmittance 75%, bar code 124.

Typical Product Data



Typical assay curve. The data was generated using a white OptiplatTM-384 microplate and the EnVision[®] Reader. The curve was obtained by mixing streptavidin acceptor and streptavidin donor beads with increasing concentrations of biotinylated mIgG. The EC₅₀ was measured from the curve portion ranging from 0 analyte to the hook point.

Please visit our website for additional information on AlphaPlex technology at www.revvy.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.

AP125TB-R Rev01

revvity