

AlphaPlex™-545

TruHits Kit

Product number: AP900TB-D **Lot Number:** 3307342

Product Format: AP900TB-D: 1 000 assay points

AP900TB-M: 10 000 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: 6/26/2024 **Document version:** 1

Kit Components

Component	AP900TB-D	AP900TB-M
AlphaPlex 545 BSA-biotin Acceptor beads stored in PBS, 0.05% Kathon, pH 7.2	1 x 50 µL at 5 mg/mL	1 x 500 µL at 5 mg/mL
Streptavidin Alpha Donor beads stored in 25 mM Hepes, 100 mM NaCl, 0.05% Kathon, pH 7.4	1 x 50 µL at 5 mg/mL	1 x 500 µL at 5 mg/mL

Product Information

Description: AlphaPlex-545 Biotin-Acceptor Beads and streptavidin donor beads at 5 mg/mL in PBS pH 7.2 supplemented with 0.05% Kathon as a preservative.

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

Storage: Store in the dark at 4°C.

Application: This product is designed for use as a tool to detect interference in AlphaPlex assays or as internal controls in Alpha assays

Quality Control

Alpha maximum signal, minimum signal are based on a one point assay performed on an EnVision® instrument. Maximum signal represents an average of the 18 wells. We certify that these results meet our requirements.

Maximum Signal: 2,441,702 counts

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

AP900TB-R Rev 01

The logo for Revvity, featuring the word "revvity" in a lowercase, sans-serif font.

Revvity, Inc.
940 Winter Street
Waltham, MA 02451 USA

(800) 762-4000 www.revivity.com

For a complete listing of our global offices, visit www.revivity.com
Copyright ©2023, Revvity, Inc. All rights reserved.