

AlphaScreen®

Protein A Acceptor beads

Product number: 6760137M Lot Number: 3285295

Product Format: 6760137M 10000 Assay Points

6760137R 50000 Assay Points

The number of assay points is based on a final bead concentration of 20 µg/mL in a 25 µL/well reaction volume.

Manufacturing date: 06/05/2024 Document version: 1

Kit Components

Component	6760137M	6760137R
Protein A Acceptor Beads at 5 mg/mL in 0.1M Tris, 0.05% Kathon, pH 8.0	1 x 1 mL (6760137)	1 x 5 mL (6760137B)

Product Information

Antibody/Protein: The Protein A acceptor beads are coated with Protein A.

Stability: This product is stable for at least 12 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.

Quality Control

Alpha maximum signal, minimum signal and EC50 are determined using a biotinylated-rIgG titration assay performed on an EnVision® instrument. We certify that these results meet our requirements.

Maximum Signal:: 524236 counts

Minimum Signal: 393 counts

EC₅₀: 0.24 nM

Typical Product Data

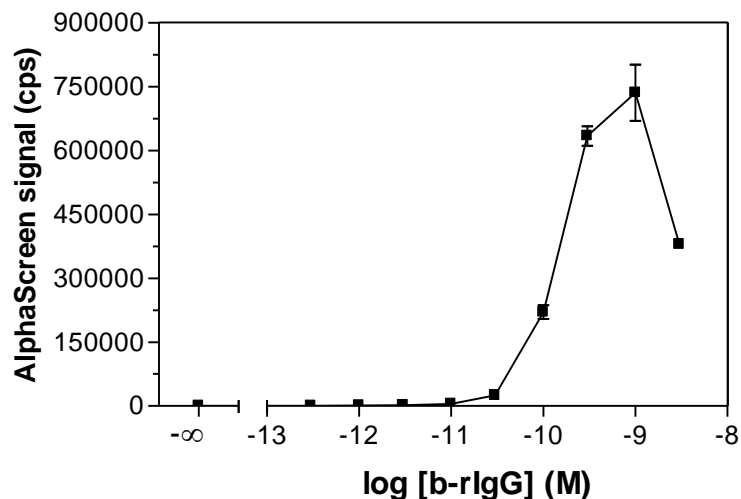


Figure 1: Biotinylated-probe titration assay 384-well biotinylated-rlgG titration curve (25 μ L final volume; Reader: Envision). Note: Alpha signal will vary depending on instrument detection protocol, incubation temperature and incubation time.

Please visit our website for additional information on AlphaLISA 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.

6760137-R Rev 01

revvity

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.