

AlphaLISA®

## Anti-Hemagglutinin (HA) Acceptor Beads

**Product number:** AL170C **Lot Number:** 3216855

---

**Material provided:** AlphaLISA Anti-HA Acceptor Beads at 5 mg/mL in PBS pH 7.2 supplemented with 0.05% Kathon as a preservative.

**Product Format:** AL170C: 250 µg, 50 µL, 500 assay points  
AL170M: 5 mg, 1 mL, 10 000 assay points  
AL170R: 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:** September 27, 2023 **Document version:** 1

---

### Product Information

**Application:** This product is designed for use in homogeneous AlphaLISA assays for the capture of HA.

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

**Stability:** This kit is stable for at least 12 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

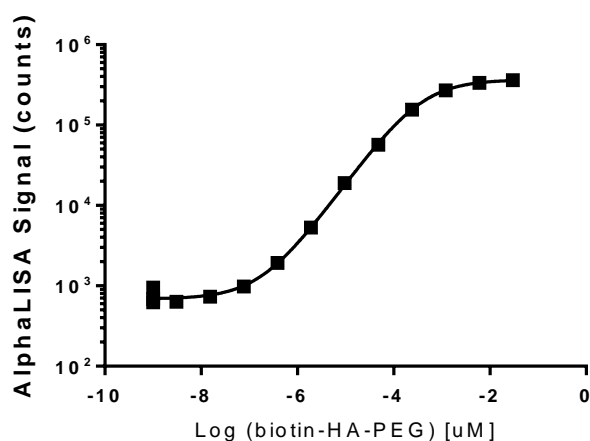
EC<sub>50</sub>: 0.05 nM  
Min counts: 248 counts  
Max counts: 363,588 counts

### Recommendations

- AlphaScreen® Donor beads are light-sensitive. All Alpha assays using the Donor beads should be performed under subdued laboratory lighting (< 100 lux). Green filters (LEE 090 filters) can be applied to light fixtures.

- Sodium azide should not be added to stock solutions or assay components. Final concentrations of sodium azide higher than 0.001 % will decrease the AlphaLISA signal.
- Small volumes may be prone to evaporation. It is recommended to cover microplates with TopSeal-A Adhesive Sealing Film to reduce evaporation during incubation. Microplates are read with the TopSeal-A Film on the plate.
- Total signal varies with temperature and incubation time. For consistent results, identical incubation times and temperature should be used for all plates.
- The AlphaLISA signal is detected with an EnVision Multilabel Reader equipped with the ALPHA option using the AlphaScreen standard settings (e.g. Total Measurement Time: 550 ms, Excitation Time: 180 ms, Mirror: D640as, Emission Filter: M570w, Center Wavelength 570 nm, Bandwidth 100 nm, Transmittance 75%).

### Typical Product Data



\* The EC50 value was determined following a non-linear regression analysis using the sigmoidal dose-response curve model with variable slope. Only assay points up to the maximum signal were used for EC50 determination.

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

AL170-R Rev01

revvity

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

(800) 762-4000 [www.revvy.com](http://www.revvy.com)

For a complete listing of our global offices, visit [www.revvy.com](http://www.revvy.com)  
Copyright ©2023, Revvity, Inc. All rights reserved.