

AlphaLISA®

Anti-Human IgE Acceptor Beads

Product number: AL176C **Lot Number:** 3292448

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

Product Format: AL176C: 250 µg, 50 µL, 500 assay points

AL176M: 5 mg, 1 mL, 10 000 assay points

AL176R: 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: April 24, 2024 **Document version:** 1

Product Information

Application: This product is intended for use in homogenous Alpha assays to capture human IgE.

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

Stability: This kit is stable for at least 6 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: 0.18 nM
Min counts: 892 counts
Max counts: 249908 counts

Titration Assay (Quality Control Procedure)

This protocol provides a means to verify product performance. The following reagents and materials are recommended.

Item	Suggested source
White OptiPlate™-384	Revvity Inc.
TopSeal™-A Plus Adhesive Sealing Film	Revvity Inc.
EnVision®-Alpha Reader	Revvity Inc.
AlphaScreen® Streptavidin coated Donor Beads	Revvity Inc.
Biotin-human IgE	Revvity Inc.
AlphaLISA Immunoassay Buffer 10X	Revvity Inc.

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.
- Spin down tubes briefly before use to improve recovery of content (2,000 x g, 10-15 sec). Resuspend all reagents by vortexing before use.
- 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%).

Protocol

- 1) Preparation of 1X AlphaLISA Immunoassay Buffer:
Add 1 mL of 10X AlphaLISA Immunoassay Buffer to 9 mL H₂O.
- 2) Preparation probe dilutions:
 - a. Reconstitute the 2.35 µg lyophilized biotin hIgE in 50 µL Milli-Q H₂O to make 250 nM stock solution.
 - a. Prepare standard dilutions as follows in 1X AlphaLISA Immunoassay Buffer (change tip between each standard dilution):

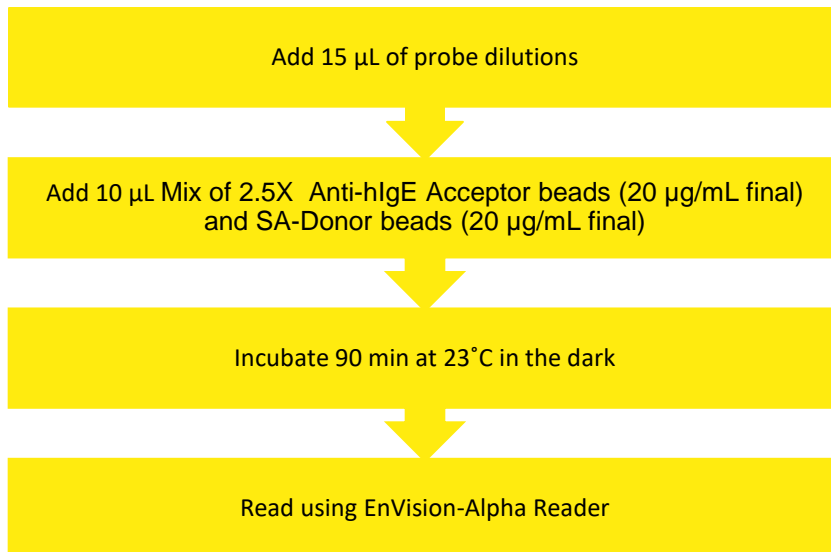
Tube	Volume of Probe	Volume of buffer (µL)	[hIgE] (M) (1.7X)	[hIgE] (pM)
A	10 µL of 250nM hIgE	140	1.7E-8	16,667
B	60 µL of tube A	140	5.0E-9	5,000
C	60 µL of tube B	120	1.7E-9	1,667

D	60 µL of tube C	140	5.0E-10	500
E	60 µL of tube D	120	1.7E-10	167
F	60 µL of tube E	140	5.0E-11	50
G	60 µL of tube F	120	1.7E-11	17
H	60 µL of tube G	140	5.0E-12	5
I	60 µL of tube H	120	1.7E-12	1.7
J	60 µL of tube I	140	5.0E-13	0.5
K	60 µL of tube J	120	1.7E-13	0.17
L	60 µL of tube K	140	5.0E-14	0.05
M	0	100	0	0
N	0	100	0	0
O	0	100	0	0
P	0	100	0	0

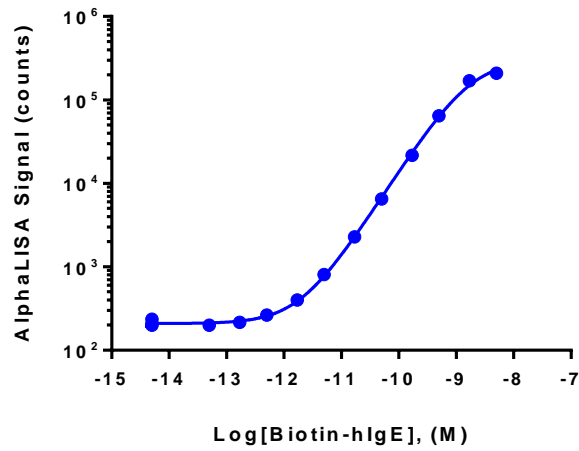
3) Preparation of the mix of 2.5X anti-hIgE Acceptor beads (50 µg/mL) and Streptavidin (SA) Donor beads (50 µg/mL) (50 µg/mL)

Add 10 µL of 5 mg/mL AlphaLISA anti-hIgE Acceptor beads and 10 µL of 5 mg/mL Streptavidin (SA) Donor beads into 980 µL of 1X AlphaLISA Immunoassay Buffer.

5) In a OptiPlate-384 microplate:



Typical Product Data



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. REVIVITY 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.

AL176-R Rev01

revvity