

AlphaScreen®

Glutathione Donor Beads

Product number: 6765301 **Lot Number:** 3299400

Material provided: Glutathione Donor Beads at 5 mg/mL in 25 mM Hepes, 0.1% Tween-20, 0.05% Kathon, pH 7.4

Product Format: 6765300: 1 mg*, packaged as 1 x 200 µL

6765301: 5 mg, packaged as 1 x 1 mL

6765302: 25 mg, packaged as 1 x 5 mL

* Note: 1mg translates into 2,000 assay points based on a final bead concentration of 20 µg/mL in a 25 µL/well reaction volume.

Manufacturing date: June 24, 2024 **Document version:** 1

Product Information

Intended Use: AlphaScreen® Glutathione Donor beads specifically bind to GST-tagged proteins. This product should be used in conjunction with an AlphaScreen® or AlphaLISA® Acceptor bead.

Stability: This product is stable for at least 24 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.

Recommended Use: 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.

Quality Control

Alpha maximum signal and background signal are determined using anti-GST acceptor beads with 100 µM GST, and performed on an EnVision® instrument. Current lot was assayed in parallel with a previous (Reference) Glutathione donor beads lot to ensure consistency. A Ratio to Reference (RTR) was calculated from these results We certify that these results meet our requirements.

<u>TEST GST</u>	<u>RESULTS</u>	<u>REFERENCE</u>	<u>RTR</u>
Maximum signal	219273	7828	28.01
Background	1854	9818	

Please visit our website for additional information on AlphaLISA 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.

676530X-R Rev 01

The Revvity logo consists of the word "revvity" in a lowercase, sans-serif font. The letters are black and have a modern, clean appearance.

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

(800) 762-4000 www.revvy.com

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