

AlphaScreen®

Biotinylated-GST (Glutathione S-Transferase)

Product number: 6760305M **Lot Number:** 3330631

Product Format: 6760305M: 10 000 assay points

The number of assay points is based on a final probe concentration of 3 nM in a 25 µL/well

Manufacturing date: 10/24/2024 **Document version:** 1

Kit Components

Component	6760305M
Biotinylated-GST at 0.5 µM in 25mM HEPES, 100mM NaCl, 0.05% Kathon, 0.05% Tween 20, pH 7.4	1 x 1.5 mL (6760305)

Product Information

Application: The biotinylated-GST probe is designed to be used with AlphaScreen® GST detection kit (6760603). Due to the length of the linker used for the biotinylation, the biotinylated-GST cannot be recognized by AlphaScreen® GSH donor beads (6765300, 6765301, 6765302 and 6765303).

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 or lower.

Quality Control

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

Maximum Signal: 814,320 counts
Minimum Signal: 798 counts
EC50: 0.63 nM

Typical Product Data

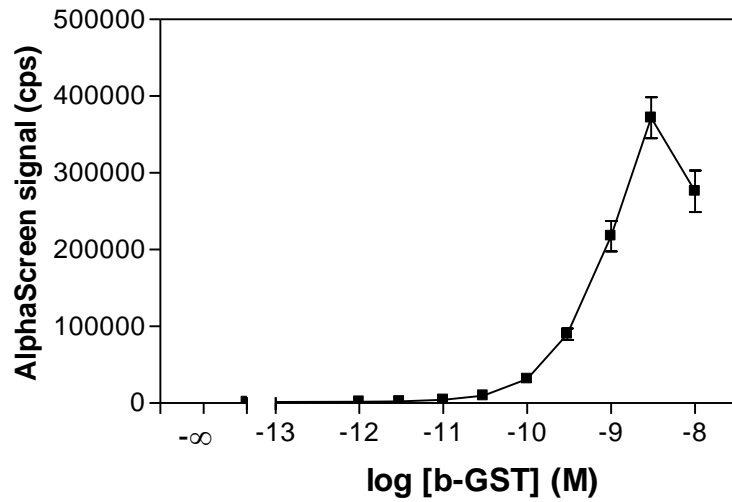


Figure 1: Biotinylated-probe titration assay 384-well biotinylated-GST titration curve (25 μ L final volume; Reader: Envision). Note: Alpha signal will vary depending on instrument detection protocol, incubation temperature and incubation time.

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

6760305-R Rev 01

revvity

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.