

AlphaScreen®

GST (Glutathione S-Transferase) Detection Kit

Product number: 6760603C

Lot Number: 3404185

Product Format: 6760603C: 500 assay points

6760603M: 10 000 assay points

6760603R: 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: May 22nd, 2025

Document version: 1

Kit Components

Component	6760603C	6760603M	6760603R
Anti-GST (Glutathione S-Transferase) Acceptor Beads at 5 mg/mL in 25 mM Hepes, 100 mM NaCl, 0.05% Kathon, pH 7.4	1 x 50 µL (6760106)	1 x 1 mL (6760107)	1 x 5 mL (6760107B)
Streptavidin Donor Beads at 5 mg/mL in 25 mM Hepes, 100 mM NaCl, 0.05% Kathon, pH 7.4	1 x 50 µL (6760001)	1 x 1 mL (6760002)	1 x 5 mL (6760002B)
Biotinylated-GST at 0.5 µM in 25mM Hepes, 100mM NaCl, 0.05% Kathon, 0.05% Tween 20, pH 7.4	1 x 50 µL (6760304)	1 x 50 µL (6760304)	1 x 50 µL (6760304)
10x Buffer : 250 mM Hepes, 1 M NaCl, 0.05% Kathon, pH 7.4	1 x 1.5 mL (6760020G)	1 x 1.5 mL (6760020G)	1 x 1.5 mL (6760020G)

Product Information

Antibody/Protein: The GST antibody is a goat polyclonal antibody that recognized Glutathione S-Transferase [Schistosoma japonicum].

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.

Storage: Store undiluted at 4°C protected from light. Freeze-thaw is not recommended and 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, 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::	405691 counts
Minimum Signal:	484 counts
EC ₅₀ :	0.52 nM

Recommended Assay Conditions

Note: This protocol provides a method to verify kit performance and is not representative of an assay. Sufficient biotinylated-probe and 10x buffer is provided to perform 3 titration curves in triplicate as described.

1x Buffer:	Add 500 µL 10x buffer to 4.5 mL Milli-Q® H ₂ O (or equivalent). Add 5 mg BSA (0.1% final concentration) and adjust pH to 7.4.
Acceptor Beads:	Add 5 µL Anti-GST (Glutathione S-Transferase) Acceptor beads to 245 µL 1x buffer.
Donor beads:	Add 5 µL Streptavidin Donor beads to 245 µL 1x buffer.
Biotinylated-probe:	From the 0.5 µM biotin-GST, prepare a ½ log dilution series (50 nM to 5 pM) in 1x buffer. Include a buffer only control.

Titration Protocol

To a white opaque OptiPlate-384:

- 1) Add 5 µL biotin-GST dilutions (from lowest to highest concentration).
- 2) Add 10 µL of 1x buffer
- 3) Add 5 µL of Anti-GST (Glutathione S-Transferase) Acceptor beads.

Incubate in the dark at room temperature for 30 minutes

- 4) Add 5 µL of Streptavidin Donor beads.

Incubate in the dark at room temperature for 60 minutes and analyze on your Alpha capable detection reader.

Typical Product Data

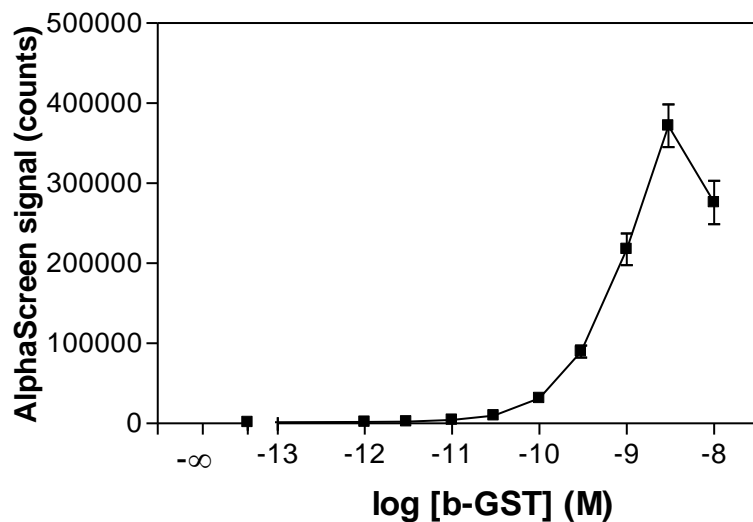


Figure 1: Biotinylated-probe titration assay 384-well biotinylated-GST titration curve (25 μ L final volume; Reader: Envision). Note: Alpha signal will vary depending on instrument detection protocol, incubation temperature and incubation time.

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.

6760603-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.