

LANCE® Ultra

ULight™ -labeled Anti-GST**Product number:** TRF0104-D **Lot Number:** 3221094**Product Format:**
TRF0104-D: 1 nmole (200µL)
TRF0104-M: 10 nmoles (2 x 1mL)
TRF0104-R: 100 nmoles (2 x 10mL)**Manufacturing date:** 10/30/2023 **Document version:** 1**Product Information****Storage Buffer:** 50 mM Tris-HCl (pH 7.4), 0.9% NaCl, 0.1% BSA and 0.05% sodium azide as preservative
Molecular Weight: 160 000
Stability: This product is stable for at least **18 months** from the manufacturing date when stored in its original packaging and the recommended storage conditions.
Storage Conditions: Store at -20°C. Store protected from light. Repeated freezing and thawing should be avoided.**Quality Control**

The QC release specifications are based on spectrophotometric analysis of the labeled antibody. We certify that results meet our quality release criteria.

Labeling Ratio: 1.7 (dye molecule/Ab)
Concentration: 800 µg/mL (5 µM)**Recommended Assay Conditions****GST tag detection assay****SUGGESTED METHOD:**
(*Specific applications might require optimization*)Reagent Preparation

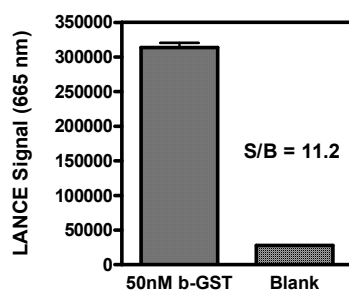
- Prepare 1X Detection Buffer: dilute 0.5 mL of 10X Detection Buffer with 4.5 mL of dH₂O.
- Prepare 2X Europium-labeled streptavidin: dilute stock solution to a concentration of 20 nM in 1X Detection Buffer.
- Prepare a 4X biotin-GST solution: dilute biotin-GST to a concentration of 200 nM in 1X Detection Buffer.
- Prepare a 4X ULight-anti GST antibody solution: dilute ULight-anti GST to a concentration of 40 nM in 1X Detection Buffer.

Protocol

- Pipet 10 µL of Europium-labeled streptavidin into a 384-well white OptiPlate™-384 (10 nM final concentration).
- Add 5 µL of biotin-GST solution (50 nM final concentration).
- Add 5 µL of *ULight*-anti GST antibody solution (10 nM final concentration).
- Cover plate with TopSeal-A and incubate 60 min at 23°C.
- Remove TopSeal-A and read in TR-FRET mode at 665 nm (excitation at 320 or 340 nm).

Typical Product Data

Detection of biotin-GST using
ULight-anti GST and
Eu (W1024)-streptavidin obtained using the EnVision® Multilabel
Reader:



Suggested Materials

- Antibody: *ULight*™ - anti GST
- Substrate: biotin-GST
- Detection: Eu--W1024-labeled streptavidin
- Detection Buffer: LANCE® Detection Buffer, 10X
- Plate: OptiPlate™-384, white
- TopSeal™-A

Supplier

Revvity Inc
Revvity Inc
Revvity Inc
Revvity Inc
Revvity Inc
Revvity Inc

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.

TRF0104-R Rev01

revvity

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

(800) 762-4000 www.revvity.com

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