

LANCE® Ultra

ULight™ -labeled Poly GAT

Product number: TRF0101-D Lot Number: 3368795

Product Format: TRF0101-D: 1 nmole (100µL)
TRF0101-M: 10 nmoles (1mL)

Manufacturing date: 01/09/2025 Document version: 1

Product Information

Phosphorylation Motif: [EAY(1:1:1)]n
Storage Buffer: 50 mM Tris-HCl (pH 7.4), 0.9% NaCl, 0.1% BSA and 0.05% sodium azide as preservative
Molecular Weight: 35 000
Stability: This product is stable for at least **10 months** from the manufacturing date when stored in its original packaging and the recommended storage conditions.
Storage Conditions: Store at -80°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: 0.6 (dye molecule/peptide)
Concentration: 350 µg/mL (10 µM)

Recommended Assay Conditions

Src kinase assay: ATP titration

SUGGESTED METHOD:

(Specific applications might require optimization)

Reagent Preparation

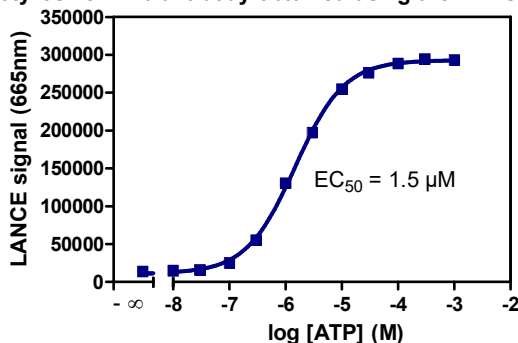
- Prepare 1X Kinase Assay Buffer: 50 mM Tris-HCl pH 7.5, 1 mM EGTA, 10 mM MgCl₂, 2 mM DTT and 0.01% Tween-20.
- Prepare a 2X c-Src enzyme solution: dilute stock enzyme to concentration of 2 nM in Kinase Assay Buffer. Keep on ice.
- Prepare a 4X ULight-poly GAT solution: dilute ULight-poly GAT to a concentration of 400 nM in Kinase Assay Buffer. Keep on ice.
- Prepare a 4X ATP solution: dilute ATP to concentrations ranging from 40 nM to 4 mM (serial half-log dilutions) in Kinase Assay Buffer. Keep on ice.
- Prepare 1X Detection Buffer: dilute 0.5 mL of 10X detection buffer with 4.5 mL of H₂O.
- Prepare a 4X Stop Solution: prepare a 40 mM EDTA solution in 1X Detection Buffer.
- Prepare a 4X Detection Mix: dilute the Eu-W1024-labeled PY20 antibody to a concentration of 8 nM in 1X Detection Buffer.

Protocol

- Pipet 5 µL of 2X c-Src enzyme into a 384-well white OptiPlate-384 (1 nM final concentration).
- Add 2.5 µL of 4X ULight-poly GAT solution (100 nM final concentration).
- Add 2.5 µL of 4X ATP solution (10 nM to 1 mM final concentrations).
- Cover plate with TopSeal-A and incubate 60 min at 23°C.
- Add 5 µL of 4X Stop Solution and incubate 5 min at 23°C.
- Add 5 µL of 4X Detection Mix (2 nM Eu-labeled PY20 antibody final concentration) and mix.
- 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

Src kinase assay using ULight-poly GAT and Eu-anti phosphotyrosine PY20 antibody obtained using the EnVision® Multilabel Reader:



Suggested Materials

- Substrate: ULight™-poly Glu, Ala, Tyr (1:1:1)
- Antibody: Eu-W1024 anti-phosphotyrosine (PY20)
- Kinase: c-Src
- Detection Buffer: LANCE® Detection Buffer, 10X
- Plate: OptiPlate™-384, white

Supplier

Revvity Inc
Revvity Inc
UpState
Revvity Inc
Revvity Inc

- TopSeal™-A

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.

TRF0101-R Rev01

The logo for Revvity, featuring the word "revvity" in a lowercase, sans-serif font.

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.