

LANCE® Ultra

ULight™-labeled CREBtide

Product number: TRF0107-M

Lot Number: 3314171

Product Format: TRF0107-D: 0.5 nmole (100µL)

TRF0107-M: 5 nmoles (1mL)

Manufacturing date: 08/14/2024

Document version: 1

Product InformationPhosphorylation Motif: RRPSYRK

Synthetic peptide derived from human cAMP Response Element Binding (CREB) Protein.
Phosphorylation site: Ser133

Storage Buffer: 50 mM Tris-HCl (pH 7.4), 0.9% NaCl, 0.1% BSA and 0.05% sodium azide as preservative

Molecular Weight: 2 740

Stability: This product is stable for at least **12 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 protein. We certify that results meet our quality release criteria.

Labeling Ratio: 1 (dye molecule/peptide)

Concentration: 13.7 µg/mL (5 µM)

Recommended Assay Conditions

MSK1: ATP titration

SUGGESTED METHOD:

*(Specific applications might require optimization)*Reagent Preparation

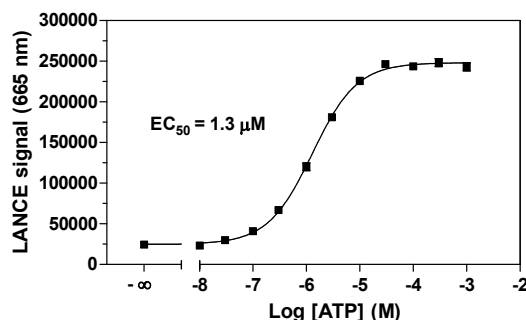
- Prepare 1X Kinase Assay Buffer: 50 mM Hepes pH 7.5, 1 mM EGTA, 10 mM MgCl₂, 2 mM DTT and 0.01% Tween-20.
- Prepare a 2X PKA solution: dilute the enzyme to a concentration of 20 pM in Kinase Assay Buffer. Keep on ice.
- Prepare a 4X mix of ULight-CREBtide: dilute ULight-CREBtide to a concentration of 200 nM in Kinase Assay Buffer.
- Prepare a 4X mix of ATP: 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-anti-phospho-CREB antibody to a concentration of 8 nM in 1X Detection Buffer.

Protocol

- Pipet 5 μ L of 2X PKA solution into a 384-well white OptiPlate™-384 (10 pM final concentration).
- Add 2.5 μ L of 4X *ULight*-CREBtide (50 nM final concentration).
- Add 2.5 μ L of 4X ATP mix (10 nM to 1 mM final concentrations).
- Cover plate with TopSeal-A and incubate 60 min at 23°C.
- Add 5 μ L of 4X Stop Solution and incubate 5 min at 23°C.
- Add 5 μ L of Detection Mix (2 nM Eu-anti-phospho-CREB 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

**PKA kinase assay using *ULight*-CREBtide
and Eu-anti-phospho-CREB antibody obtained using the EnVision® Multilabel
Reader:**



Suggested Materials

- Substrate: *ULight*™- CREBtide peptide
- Antibody: Eu- anti-phospho- CREB (Ser133)
- Kinase: PKA, catalytic subunit, recombinant
- Detection Buffer: LANCE® Detection Buffer, 10X
- Plate: OptiPlate™-384, white
- TopSeal™-A

Supplier

Revvity Inc
Revvity Inc
Upstate
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.

TRF0107-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.