

LANCE® Ultra

Eu-W1024 labeled Anti-phospho-CREB (Ser133) Antibody

Product number: TRF0200-C Lot Number: 3303443

Product Format: TRF0200-C: 1.7 µg

Manufacturing date: March 6, 2024 Document version: 1

Product Information

Antibody: Europium-labeled rabbit monoclonal antibody recognizing phospho-Ser133 of the cAMP responsive element binding (CREB) protein.

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 24 months from the manufacturing date when stored in its original packaging and at the recommended storage conditions.

Storage Conditions: Store at 4°C

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: 6.28
Concentration: 100 µg/mL

Recommended Assay Conditions

PROTEIN KINASE A: ATP TITRATION

SUGGESTED METHOD:
(*Specific applications might require optimization*)

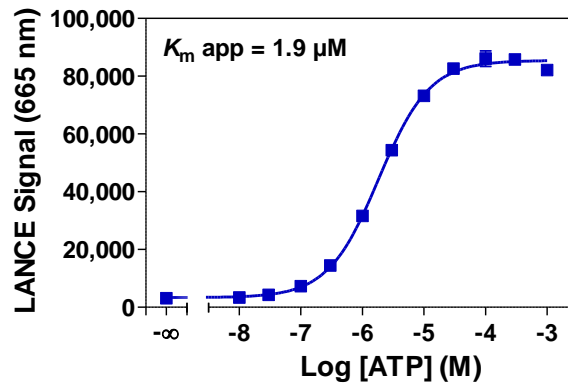
Reagent Preparation

- 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 5 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 24 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 (2.5 pM final concentration).
- Add 2.5 μ L of 4X *ULight*- CREBtide (50 nM final concentration).
- Add 2.5 μ L of 4X ATP solution (10 nM to 1 mM final concentrations).
- Cover plate with TopSeal-A and incubate 45 min at 23°C.
- Add 5 μ L of 4X Stop solution and incubate 5 min at 23°C.
- Add 5 μ L of 4X Detection Mix (2 nM Eu-anti-phospho-CREB antibody final conc.) 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



Suggested Materials

- | | |
|--|-----------------|
| • Substrate: <i>ULight</i> [™] - CREBtide) Peptide | Supplier |
| • Antibody: Eu- phospho-CREB (Ser133) Antibody | Revvity Inc |
| • Kinase: PKA, catalytic subunit, recombinant | Revvity Inc |
| • Detection Buffer: LANCE [®] Detection Buffer, 10X | Millipore |
| • Plate: OptiPlate [™] -384, white | 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.

TRF0200-C-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.