

LANCE® Ultra

Eu-W1024 labeled Anti-phospho-MEK1/MEK2 Antibody**Product number:** TRF0213-D **Lot Number:** 3341752**Product Format:** TRF0213-D: 10 µg

TRF0213-M: 100 µg

Manufacturing date: 9/19/2024 **Document version:** 1**Product Information****Antibody:** Europium-labeled rabbit polyclonal antibody recognizing phospho-Ser217/221 in human MEK1/2**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 15 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: 6.9 / 1**Concentration:** 0.625 µM, 100 µg/mL**Recommended Assay Conditions**

Detection of a Phospho-MEK1/2 Biotinylated Peptide

SUGGESTED METHOD:

(Specific applications might require optimization)

Reagent Preparation

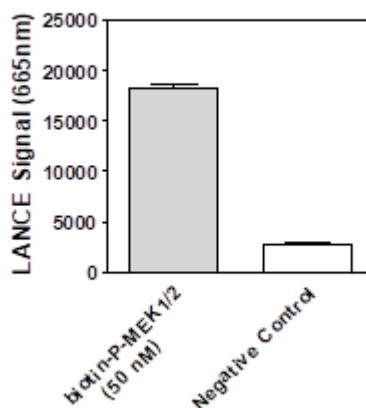
- Prepare 1X Kinase Assay Buffer: 50 mM HEPES pH 7.5, 1 mM EGTA, 10 mM MgCl₂, 2 mM DTT and 0.01% Tween-20.
- Prepare a 1X biotin-phospho-MEK1/2 (Ser217/221) Peptide solution: dilute biotin-phospho-MEK1/2 (Ser217/221) Peptide to a concentration of 50 nM in 1X Kinase Assay Buffer.
- Prepare 1X Detection Buffer: dilute 0.5 mL of 10X Detection Buffer with 4.5 mL of H₂O.
- 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 Europium-anti-phospho-MEK1/2 (Ser217/221) Antibody solution: dilute Europium-anti-phospho-MEK1/2 (Ser217/221) Antibody to a concentration of 8 nM in 1X Detection Buffer.
- Prepare a 4X *ULight*-Streptavidin solution: dilute *ULight*-Streptavidin to a concentration of 200 nM in 1X Detection Buffer.

Protocol

- Pipet 10 µL of 1X biotin-phospho-MEK1/2 (Ser217/221) Peptide solution into a 384-well white OptiPlate-384 (50 nM final concentration).
- Add 5 µL of 4X Europium-anti-phospho-MEK1/2 (Ser217/221) Antibody solution to all wells (2 nM final concentration).
- Add 5 µL of 4X *ULight*-Streptavidin (50 nM final concentration). Add buffer only to control wells
- 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 a biotinylated phospho-MEK1/2 peptide
obtained using the EnVision® Multilabel Reader:



Suggested Materials

- Detection: *ULight*[™]-Streptavidin
- Antibody: Eu- anti-phospho-MEK1/2 (Ser217/221)
- Phospho-MEK1/2 (Ser217/221) Biotinyl. Pept.
- Detection Buffer: LANCE[®] Detection Buffer, 10X
- Plate: OptiPlate[™]-384, white
- TopSeal[™]-A

Supplier

Revvity Inc
Revvity Inc
CST
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.

TRF0213-R Rev01

revvity