

HTRF® Research Reagents

HTRF MAb Anti p-ELK-1-Eu - 500 pts
Product number: 61P11KAE

Lot number: 05RA

Product characteristics:

| | |
|---|----------------------------|
| • Packaging Unit | 500 tests/vial |
| • Conditioning buffer | Hepes 0.05M pH7 + 0.1% BSA |
| • Appearance | Frozen form |
| • Storage conditions | -20 °C |
| • Expiration date | 2026.01.18 |
| • Volume per vial | 250.00 µL |
| • Active Moiety Concentration (Ab) | 3.00 µg/mL |

Weight indications:

| | Quantity per vial | MW (Da) |
|--------------------|-------------------|---------|
| Active Moiety (Ab) | 0.75 µg | 150 000 |

Recommended quantity per well*:

| | in 384 well low volume plate (20µL) |
|----------------------|-------------------------------------|
| Active Moiety (Ab) | 1.50 ng |
| Number of tests/vial | 500 |

*Actual amount per well should be optimized according to assay conditions

Please note

KF can play an essential role in the lanthanide cryptate protection by preventing the action of possible quenchers contained in the assay. KF is mandatory for HTRF assays using Europium cryptate. KF is generally used at a final concentration of 100 to 400 mM, and is added to the conjugate working solutions or dispensed in a separate step, just before the readout. The expiration of batch 05RA has been extended to 2026-01-18 after validation by QC

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.