

phosph MEK1/2-K  
Product reference :

PAb Anti-phospho MEK1/2-Eu cryptate - 10,000 tests  
61P17KAZ

**Product characteristics :**

- Lot Number \_\_\_\_\_ 26B
- Packaging Unit \_\_\_\_\_ 10 000 tests/vial
- Conditioning buffer \_\_\_\_\_ P04 0.1M pH7 + 0.1% BSA
- Appearance \_\_\_\_\_ Frozen form
- Storage conditions \_\_\_\_\_ <-16 °C
- Expiration date \_\_\_\_\_ 2028.05.20
- Volume per vial \_\_\_\_\_ 0.25 mL
- Active Moiety Concentration (Ab) \_\_\_\_\_ 120 µg/mL

**Weight indications :**

|                    | Quantity per vial | MW (Da) |
|--------------------|-------------------|---------|
| Active Moiety (Ab) | 30.00 µg          | 150 000 |

**Recommended quantity per well\* :**

|                      | in 384 well<br>low volume<br>plate (20µL) |
|----------------------|---|
| Active Moiety (Ab)   | 3.00 ng                                   |
| Number of tests/vial | 10 000                                    |

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

**PLEASE NOTE**

Lot to lot variability of the HTRF conjugate may be observed according to the antiphospho MEK1/2 polyclonal antibody source whose batch to batch phosphoreactivity may change .

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.

**Further information as well as a specific application note regarding conjugate-related calculations are available on our Life Sciences web site: [www.cisbio.com](http://www.cisbio.com)**