

phosph P44/42-K
Product reference :

MAb Anti-phosphop44/42-Eu cryptate - 10,000 tests
61P10KAZ

Product characteristics :

- Lot Number _____ 008
- Packaging Unit _____ 10 000 tests/vial
- Conditioning buffer _____ Hepes 50 mM pH7 + 0.1% BSA
- Appearance _____ Frozen form
- Storage conditions _____ -20 °C
- Expiration date _____ 2023.01.19
- Volume per vial _____ 250 µL
- Active Moiety Concentration (Ab) _____ 84 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Ab)	21.00 µg	150 000

Recommended quantity per well* :

	in 384 well low volume plate (20µL)
Active Moiety (Ab)	2.10 ng
Number of tests/vial	10 000

*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 date of batch 08 has been extended to 2023-01-19 after validation by QC

Further information as well as a specific application note regarding conjugate-related calculations are available on our Life Sciences web site: www.cisbio.com