

phosph PKC-K
Product reference :

MAb Anti-phospho PKC substrate-Eu cryptate - 10,000 tests
61P03KAZ

Product characteristics :

- Lot Number_____ 07A
- Packaging Unit_____ 10 000 tests/vial
- Conditioning buffer_____ hepes 50 mM pH7 + 0.1% BSA
- Appearance_____ Frozen form
- Storage conditions_____ <-16 °C
- Expiration date_____ 2028.07.31
- Volume per vial_____ 250 µL
- Active Moiety Concentration (Ab)_____ 116 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Ab)	29.00 µg	150 000

Recommended quantity per well* :

	in 384 well low volume plate (20µL)
Active Moiety (Ab)	2.90 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.

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