

phosph MEK1/2-K
Product reference :

PAb Anti-phospho MEK1/2-Eu cryptate - 100,000 tests
61P17KAY

Product characteristics :

- Lot Number	28A
- Packaging Unit	100 000 tests/vial
- Conditioning buffer	P04 0.1M Ph7 + 0.1% BSA
- Appearance	Frozen form
- Storage conditions	<-16 °C
- Expiration date	2029.07.21
- Volume per vial	2.5 mL
- Active Moiety Concentration (Ab)	120 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Ab)	0.30 mg	150 000

Recommended quantity per well* :

	in 384 well low volume plate (20µL)
Active Moiety (Ab)	3.00 ng
Number of tests/vial	100 000

*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