

phosph p53-K
Product reference :

MAb Anti-phospho p53-K - 10,000 tests
61P08KAZ

Product characteristics :

- Lot Number _____ 008
- Packaging Unit _____ 10 000 tests/vial
- Conditioning buffer _____ Hepes buffer 0.05M pH7+ 0.1% BSA
- Appearance _____ Frozen form
- Storage conditions _____ <-16 °C
- Expiration date _____ 2026.11.16
- Volume per vial _____ 2.5 mL
- Active Moiety Concentration (Ab) _____ 100 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Ab)	25.00 µg	150 000

Recommended quantity per well* :

	in 384 well low volume plate (20µL)
Active Moiety (Ab)	2.50 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