

**phosph p53-K**  
Product reference :

MAb Anti-phospho p53-Eu cryptate - 500 tests  
**61P08KAE**

**Product characteristics :**

- Lot Number \_\_\_\_\_ 08C
- Packaging Unit \_\_\_\_\_ 500 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 \_\_\_\_\_ 0.25 mL
- Active Moiety Concentration (Ab) \_\_\_\_\_ 5 µg/mL

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (Ab)	1.25 µg	150 000

**Recommended quantity per well\* :**

	in 384 well low volume plate (20µL)
Active Moiety (Ab)	2.50 ng
Number of tests/vial	500

\*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](http://www.cisbio.com)**