

phosph p53-K
Product reference :

anti phospho P53 cryptate conjugate
61P08KAZ

Product characteristics :

- Lot Number_____ 06
- Packaging Unit_____ 10 000 tests/vial
- Conditioning buffer_____ Phosphate 0.1M pH7 + 0.1% BSA + 0.1% Tween
- Appearance_____ Frozen form
- Storage conditions_____ <-16 °C
- Expiration date_____ 2024.03.25
- Volume per vial_____ 330 µL
- Active Moiety Concentration (Ab)_____ 76.5 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 HTRF web site: www.htrf.com/products