

Batch information

phosph p53-K

HTRF MAb Anti p-p53-Eu - 500 pts

Product reference :

61P08KAE

Product characteristics :

- Lot Number	10A
- Packaging Unit	500 tests/vial
- Conditioning buffer	Hepes buffer 0.05M pH7+ 0.1% BSA
- Appearance	Frozen form
- Storage conditions	<-16 °C
- Expiration date	2032.01.05
- Volume per vial	0.25 mL
- Active Moiety Concentration (Ab)	5.00 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Ab)	1 250.00 ng	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

