

ab h3k27me1-K
Product reference :

Mab Anti histone H3K27 Me1-Eu(K) - 500 tests
61KC1KAE

Product characteristics :

- Lot Number _____ 04A
 - Packaging Unit _____ 500 tests/vial
 - Conditioning buffer _____ Hepes 50mM pH7 0.1% BSA
 - Appearance _____ Frozen form
 - Storage conditions _____ -20 °C
 - Expiration date _____ 2029.02.17
 - Volume per vial _____ 50 µL
 - Active Moiety Concentration (ab) _____ 22.5 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (ab)	1.13 µg	150 000

Recommended quantity per well* :

	in 384 well low volume plate (20µL)
Active Moiety (ab)	2.25 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