

ab h3k27me1-K
Product reference :Mab Anti histone H3K27 Me1-Eu(K) - 10,000 tests
61KC1KAD**Product characteristics :**

- Lot Number	04A
- Packaging Unit	10 000 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	1000 µL
- Active Moiety Concentration (ab)	22.5 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (ab)	22.50 µg	150 000

Recommended quantity per well* :

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