

ab h3k27me2-K
Product reference :

Mab Anti histone H3K27 Me2-Eu(K) - 500 tests
61KC2KAE

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 _____ 2026.09.20
- 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

Further information as well as a specific application note regarding conjugate-related calculations are available on our Life Sciences web site: www.cisbio.com