

**ab h3k27me2-K**  
Product reference :

Mab Anti histone H3K27 Me2-Eu(K) - 10,000 tests  
**61KC2KAD**

**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 \_\_\_\_\_ 2026.09.20  
- Volume per vial \_\_\_\_\_ 1 mL  
- 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

**Further information as well as a specific application note regarding conjugate-related calculations are available on our Life Sciences web site: [www.cisbio.com](http://www.cisbio.com)**