

**ab h3k9me1-K**  
Product reference :

Mab Anti histone H3K9 Me1-Eu(K) - 500 tests  
**61KB1KAE**

**Product characteristics :**

- Lot Number\_\_\_\_\_ 04A  
- Conditioning buffer\_\_\_\_\_ HEPES 50mM pH7 +0.1% BSA +  
Tween 20 0.1%  
- Appearance\_\_\_\_\_ Frozen form  
- Storage conditions\_\_\_\_\_ -20 °C  
- Expiration date\_\_\_\_\_ 2027.05.09  
- 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

**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](http://www.cisbio.com)***