

ab h3k9me0-K
Product reference :

Mab Anti histone H3K9 Me0-Eu(K) - 10,000 tests
61KB0KAD

Product characteristics :

- Lot Number _____ 02A
- Conditioning buffer _____ HEPES 50mM pH7 + 0.1% BSA
- Appearance _____ Frozen form
- Storage conditions _____ -20 °C
- Expiration date _____ 2030.01.04
- Volume per vial _____ 1 mL
- Active Moiety Concentration (ab) _____ 15 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (ab)	15.00 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