

ab h3k36me2-K
Product reference :

Mab Anti histone H3K36 Me2-Eu(K) - 500 tests
61KD2KAE

Product characteristics :

- Lot Number_____ 05B
- Conditioning buffer_____ Hepes 50mM pH7 + 0.1% BSA
- Appearance_____ Frozen form
- Storage conditions_____ -20 °C
- Expiration date_____ 2023.05.05
- 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 HTRF web site: www.htrf.com/products