

ab h3k36me2-K
Product reference :Mab Anti histone H3K36 Me2-Eu(K) - 1mg
61KD2KAH**Product characteristics :**

- Lot Number	004B
- Packaging Unit	4 444 444 tests/vial
- Conditioning buffer	Hepes 50mM pH7 + 0.1% BSA
- Appearance	Frozen form
- Storage conditions	-20 °C
- Expiration date	2023.06.20
- Volume per vial	10 mL
- Active Moiety Concentration (ab)	100 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (ab)	1.00 mg	150 000

Recommended quantity per well* :

	in 384 well low volume plate (20µL)
Active Moiety (ab)	2.25 ng
Number of tests/vial	4 444 444

*Actual amount per well should be optimized according to assay conditions

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