

ab h3k36me1-K
Product reference :

Mab Anti histone H3K36 Me1-Eu(K) - 10,000 tests
61KD1KAD

Product characteristics :

- Lot Number	05A
- Conditioning buffer	Phosphate 0.1M pH7 + 0.1 % BSA
- Appearance	Frozen form
- Storage conditions	-20 °C
- Expiration date	2026.03.09
- 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

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

