

MAB Anti 6HIS-Eu cryptate Gold MAb Anti-6HIS-Eu cryptate Gold - 1 mg - CUSTOM LABELING
Product reference : **61HI2KAH**

Product characteristics :

- Lot Number_____ 07A
- Conditioning buffer_____ P04 100 mM pH7 + stabilizers +
BSA 0.5%
- Appearance_____ Frozen form
- Storage conditions_____ <-16 °C
- Expiration date_____ 2029.02.11
- Volume per vial_____ 5.295 mL
- Active Moiety Concentration (HIS)_____ 190 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (HIS)	1.00 mg	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