

Batch information

MAb Anti 6HIS-Eu cryptate Gold

MAb Anti-6HIS-Eu cryptate Gold - 20,000 tests

Product reference :

61HI2KLB

Product characteristics :

- Lot Number	16A
- Packaging Unit	20 000 tests/vial
- Conditioning buffer	P04 100 mM pH7 + stabilizers + BSA 0.5%
- Appearance	Lyophilized form
- Storage conditions	2-8 °C
- Expiration date	2024.06.09
- Reconstitute each vial with	1 mL distilled H ₂ O
- Active Moiety Concentration (HIS)	30.00 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (HIS)	30.00 µg	150 000

Recommended quantity per well* :

	in 384 well low volume plate (20µL)
Active Moiety (HIS)	1.50 ng
Number of tests/vial	20 000

*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

