

MAB Anti 6His Tb cryptate Gold MAB Anti-6HIS Tb cryptate Gold - 5,000 tests
 Product reference : **61HI2TLA**

Product characteristics :

- Lot Number _____ 10A
- Packaging Unit _____ 5 000 tests/vial
- Conditioning buffer _____ P04 100 mM pH7 + stabilizers +
BSA 0.5%
- Appearance _____ Lyophilized form
- Storage conditions _____ 2-8 °C
- Expiration date _____ 2022.01.22
- Reconstitute each vial with _____ 0.25 mL distilled H2O
- Active Moiety Concentration (HIS) _____ 21 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (HIS)	5.25 µg	150 000

Recommended quantity per well* :

	in 384 well low volume plate (20µL)
Active Moiety (HIS)	1.05 ng
Number of tests/vial	5 000

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

PLEASE NOTE

The use of KF is not compulsory in HTRF assays using Lumi4-Tb conjugate.

Further information as well as a specific application note regarding conjugate-related calculations are available on our Life Sciences web site: www.cisbio.com