

MAB Anti 6His Tb cryptate Gold MAb Anti-6HIS Tb cryptate Gold - 1 mg - CUSTOM LABELING
 Product reference : **61HI2TAH**

Product characteristics :

- Lot Number _____ 10A
- Packaging Unit _____ 952 380 tests/vial
- Conditioning buffer _____ P04 100 mM pH7 + BSA 0.1%
- Appearance _____ Frozen form
- Storage conditions _____ <-16 °C
- Expiration date _____ 2029.01.20
- Volume per vial _____ 4 mL
- Active Moiety Concentration (HIS) _____ 250 µg/mL

Weight indications :

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

Recommended quantity per well* :

	in 384 well low volume plate (20µL)
Active Moiety (HIS)	1.05 ng
Number of tests/vial	952 380

*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