

**MAB Anti 6His Tb cryptate Gold**    MAb Anti 6HIS-Tb cryptate Gold - 1,000 tests  
 Product reference :                      **61HI2TLF**

**Product characteristics :**

- Lot Number\_\_\_\_\_ 09A
- Packaging Unit\_\_\_\_\_ 1 000 tests/vial
- Conditioning buffer\_\_\_\_\_ P04 100 mM pH7 + stabilizers +  
BSA 0.5%
- Appearance\_\_\_\_\_ Lyophilized form
- Storage conditions\_\_\_\_\_ 2-8 °C
- Expiration date\_\_\_\_\_ 2022.09.01
- Reconstitute each vial with\_\_\_\_\_ 0.25 mL distilled H2O
- Active Moiety Concentration (HIS)\_\_\_\_\_ 4.2 µg/mL

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (HIS)	1.05 µg	150 000

**Recommended quantity per well\* :**

	in 384 well low volume plate (20µL)
Active Moiety (HIS)	1.05 ng
Number of tests/vial	1 000

\*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](http://www.cisbio.com)**

