

**MAB Anti 6His Tb cryptate Gold**    MAb Anti-6HIS Tb cryptate Gold - 20,000 tests  
Product reference :                      **61HI2TLB**

**Product characteristics :**

- Lot Number \_\_\_\_\_ 10A
- 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 \_\_\_\_\_ 2022.01.22
- Reconstitute each vial with \_\_\_\_\_ 1 mL distilled H2O
- Active Moiety Concentration (HIS) \_\_\_\_\_ 21 µg/mL

**Weight indications :**

|                     | Quantity per vial | MW (Da) |
|---------------------|-------------------|---------|
| Active Moiety (HIS) | 21.00 µg          | 150 000 |

**Recommended quantity per well\* :**

|                      | in 384 well<br>low volume<br>plate (20µL) |
|----------------------|---|
| Active Moiety (HIS)  | 1.05 ng                                   |
| Number of tests/vial | 20 000                                    |

\*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)**