

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

**Product characteristics :**

- Lot Number	08A
- Packaging Unit	952 380 tests/vial
- Conditioning buffer	P04 100 mM pH7 + 0.1% BSA
- Appearance	Frozen form
- Storage conditions	<16 °C
- Expiration date	2027.09.25
- Volume per vial	2.36 mL
- Active Moiety Concentration (HIS)	425 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](http://www.cisbio.com)**