

## Batch information

**MAb Anti 6His Tb cryptate Gold**

MAb Anti-6HIS-Tb cryptate Gold - 1,000 tests

Product reference :

**61HI2TLF**

### Product characteristics :

- Lot Number	17A
- 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	2024.03.31
- Reconstitute each vial with	0.25 mL distilled H2O
- Active Moiety Concentration (HIS)	4.00 µg/mL

### Weight indications :

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

### Recommended quantity per well\* :

	in 384 well low volume plate (20µL)
Active Moiety (HIS)	1.00 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)**

