

**MAB 6HIS-Tb**

MAB Anti 6HIS-Tb cryptate - 1 mg - CUSTOM LABELING  
**61HISTAH**

Product reference :

**Product characteristics :**

- Lot Number\_\_\_\_\_ 07A
- Packaging Unit\_\_\_\_\_ 555 555 tests/vial
- Conditioning buffer\_\_\_\_\_ Phosphate 0.1M pH7 + 0.1% BSA +  
0.1% Tween
- Appearance\_\_\_\_\_ Frozen form
- Storage conditions\_\_\_\_\_ - 16 °C
- Expiration date\_\_\_\_\_ 2025.04.27
- Volume per vial\_\_\_\_\_ 4 mL
- Active Moiety Concentration (MAB)\_\_\_\_\_ 250 µg/mL

**Weight indications :**

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

**Recommended quantity per well\* :**

	in 384 well low volume plate (20µL)
Active Moiety (MAB)	1.80 ng
Number of tests/vial	555 555

\*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.

The performance of assays using 6HIS-tag can be enhanced by incubating those at 4-8°C.

**Further information as well as a specific application note regarding conjugate-related calculations are available on our HTRF web site: [www.htrf.com/products](http://www.htrf.com/products)**