

**MAb anti HA-Tb**  
Product reference :

MAb Anti HA-Tb cryptate - 1,000 tests  
**610HATAF**

**Product characteristics :**

- Lot Number \_\_\_\_\_ 09B  
- Packaging Unit \_\_\_\_\_ 1 000 tests/vial  
- Conditioning buffer \_\_\_\_\_ Phosphate 0.1M pH7 + 0.1% BSA  
- Appearance \_\_\_\_\_ Frozen form  
- Storage conditions \_\_\_\_\_ <-16 °C  
- Expiration date \_\_\_\_\_ 2026.12.20  
- Volume per vial \_\_\_\_\_ 0.05 mL  
- Active Moiety Concentration (Mab) \_\_\_\_\_ 24 µg/mL

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (Mab)	1.20 µg	150 000

**Recommended quantity per well\* :**

	in 384 well low volume plate (20µL)
Active Moiety (Mab)	1.20 ng
Number of tests/vial	1 000

\*Actual amount per well should be optimized according to assay conditions

**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)**