

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

Monoclonal Anti HA Antibody labeled with Lumi4 Tb  
**610HATAB**

**Product characteristics :**

- Lot Number\_\_\_\_\_ 07A  
 - Packaging Unit\_\_\_\_\_ 20 000 tests/vial  
 - Conditioning buffer\_\_\_\_\_ Phosphate 0.1M pH7 + 0.1% BSA  
 - Appearance\_\_\_\_\_ Frozen form  
 - Storage conditions\_\_\_\_\_ <-16 °C  
 - Expiration date\_\_\_\_\_ 2024.09.30  
 - Volume per vial\_\_\_\_\_ 1 mL  
 - Active Moiety Concentration (Mab)\_\_\_\_\_ 24 µg/mL

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (Mab)	24.00 µg	150 000

**Recommended quantity per well\* :**

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

\*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 HTRF web site: [www.htrf.com/products](http://www.htrf.com/products)**