

MAb anti HA-Tb
Product reference :

MAb Anti HA-Tb cryptate - 5,000 tests
610HATAA

Product characteristics :

- Lot Number _____ 08A
 - Packaging Unit _____ 5 000 tests/vial
 - Conditioning buffer _____ Phosphate 0.1M pH7 + 0.1% BSA
 - Appearance _____ Frozen form
 - Storage conditions _____ <-16 °C
 - Expiration date _____ 2026.02.21
 - Volume per vial _____ 0.25 mL
 - Active Moiety Concentration (Mab) _____ 24 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Mab)	6.00 µg	150 000

Recommended quantity per well* :

	in 384 well low volume plate (20µL)
Active Moiety (Mab)	1.20 ng
Number of tests/vial	5 000

*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