

MAb anti HA-Tb
Product reference :

MAb Anti HA-Tb cryptate - 20,000 tests
610HATAB

Product characteristics :

- Lot Number _____ 08A
- Packaging Unit _____ 20 000 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 _____ 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