

MAb anti HA-Tb
Product reference :MAb Anti HA-Tb cryptate - 5,000 tests - CUSTOM LABELING
610HATAA**Product characteristics :**

- Lot Number	10A
- Packaging Unit	5 000 tests/vial
- Conditioning buffer	Phosphate 0.1M pH7 + 0.1% BSA
- Appearance	Frozen form
- Storage conditions	<-16 °C
- Expiration date	2028.09.02
- 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 Life Sciences web site: www.cisbio.com