

Anti-c-myc-Tb

MAb Anti cmyc-Tb cryptate - 20,000 tests

Product reference :

61MYCTAB**Product characteristics :**

- Lot Number	05A
- Packaging Unit	20 000 tests/vial
- Conditioning buffer	Phosphate 0.1M pH7 + 0.1 % BSA
- Appearance	Frozen form
- Storage conditions	<-16 °C
- Expiration date	2029.06.21
- Volume per vial	1 mL
- Active Moiety Concentration (MAb)	30 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (MAb)	30.00 µg	150 000

Recommended quantity per well* :

	in 384 well low volume plate (20µL)
Active Moiety (MAb)	1.50 ng
Number of tests/vial	20 000

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

PLEASE NOTE

The use of KF is not compulsory in HTRF assays using Lumi4-Tb conjugate.

Further information as well as a specific application note regarding conjugate-related calculations are available on our Life Sciences web site: www.cisbio.com