

Anti-c-myc-Tb

MAb Anti cmyc-Tb cryptate - 5,000 tests

Product reference :

61MYCTAA

Product characteristics :

- Lot Number	03B
- Packaging Unit	5 000 tests/vial
- Conditioning buffer	Phosphate 0.1M pH7 + 0.1 % BSA
- Appearance	Frozen form
- Storage conditions	-20 °C
- Expiration date	2024.05.04
- Volume per vial	0.25 mL
- Active Moiety Concentration (MAb)	30 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (MAb)	7.50 µg	150 000

Recommended quantity per well* :

	in 384 well low volume plate (20µL)
Active Moiety (MAb)	1.50 ng
Number of tests/vial	5 000

*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 HTRF web site: www.htrf.com/products