

Anti-c-myc-d2

Product reference :

MAb Anti cmyc-d2 - 5,000 tests

61MYCDAA

Product characteristics :

- Lot Number	15A
- Packaging Unit	5 000 tests/vial
- Conditioning buffer	Phosphate 0.1M pH7 + 0.1% BSA
- Appearance	Frozen form
- Storage conditions	< -16 °C
- Expiration date	2028.02.04
- Volume per vial	400 µL
- Active Moiety Concentration (MAb)	250 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (MAb)	100.00 µg	150 000

Recommended quantity per well* :

	in 384 well low volume plate (20µL)
Active Moiety (MAb)	20.00 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