

**Anti-c-myc-d2**

MAb Anti cmyc-d2 - 20,000 tests

Product reference :

**61MYCDAB**

**Product characteristics :**

- Lot Number_____	15A
- Packaging Unit_____	20 000 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_____	1.6 mL
- Active Moiety Concentration (MAb)_____	250 µg/mL

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (MAb)	0.40 mg	150 000

**Recommended quantity per well\* :**

	in 384 well low volume plate (20µL)
Active Moiety (MAb)	20.00 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 Life Sciences web site: [www.cisbio.com](http://www.cisbio.com)**