

Anti-c-myc-d2
Product reference :

Monoclonal Anti-c-myc Antibody labeled with d2
61MYCDAB

Product characteristics :

- Lot Number_____ 08A
- Packaging Unit_____ 20 000 tests/vial
- Conditioning buffer_____ Phosphate 0.1M pH7 + 0.1% BSA
- Appearance_____ Frozen form
- Storage conditions_____ <-16 °C
- Expiration date_____ 2024.06.08
- 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 HTRF web site: www.htrf.com/products