

**Anti-c-myc-d2**  
Product reference :

MAb Anti cmyc-d2 - 20,000 tests  
**61MYCDAB**

**Product characteristics :**

- Lot Number \_\_\_\_\_ 12A  
 - Packaging Unit \_\_\_\_\_ 20 000 tests/vial  
 - Conditioning buffer \_\_\_\_\_ Phosphate 0.1M pH7 + 0.1% BSA  
 - Appearance \_\_\_\_\_ Frozen form  
 - Storage conditions \_\_\_\_\_ < -16 °C  
 - Expiration date \_\_\_\_\_ 2025.05.31  
 - 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](http://www.htrf.com/products)**