

**Anti HA - d2**

MAb Anti HA-d2 - 20,000 tests

Product reference :

**610HADAB**

**Product characteristics :**

- Lot Number \_\_\_\_\_ 06B  
 - Packaging Unit \_\_\_\_\_ 20 000 tests/vial  
 - Conditioning buffer \_\_\_\_\_ Phosphate 0.1M pH7 + 0.1 % BSA  
 - Appearance \_\_\_\_\_ Frozen form  
 - Storage conditions \_\_\_\_\_ <-16 °C  
 - Expiration date \_\_\_\_\_ 2027.06.19  
 - Volume per vial \_\_\_\_\_ 1.6 mL  
 - Active Moiety Concentration (Mab) \_\_\_\_\_ 0.25 mg/mL

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (Mab)	400.00 µg	150 000

**Recommended quantity per well\* :**

	in 384 well low volume plate (20µL)
Active Moiety (Mab)	20.00 ng
Number of tests/vial	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)**