

**Anti HA - d2**

MAb Anti HA-d2 - 1,000 tests

Product reference :

**610HADAF**

**Product characteristics :**

- Lot Number\_\_\_\_\_ 06B  
- Packaging Unit\_\_\_\_\_ 1 000 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\_\_\_\_\_ 0.08 mL  
- Active Moiety Concentration (Mab)\_\_\_\_\_ 0.25 mg/mL

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (Mab)	20.00 µg	150 000

**Recommended quantity per well\* :**

	in 384 well low volume plate (20µL)
Active Moiety (Mab)	20.00 ng
Number of tests/vial	1 000

\*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)**