

Anti HA - d2

MAb Anti HA-d2 - 20,000 tests

Product reference :

610HADAB

Product characteristics :

- Lot Number _____ 07A
 - Packaging Unit _____ 20 000 tests/vial
 - Conditioning buffer _____ Phosphate 0.1M pH7 + 0.1 % BSA
 - Appearance _____ Frozen form
 - Storage conditions _____ <-16 °C
 - Expiration date _____ 2027.12.10
 - Volume per vial _____ 1.6 mL
 - Active Moiety Concentration (Mab) _____ 250 µg/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