

Anti HA - d2

MAb Anti HA-d2 - 20,000 tests

Product reference :

610HADAB**Product characteristics :**

- Lot Number	09A
- Packaging Unit	20 000 tests/vial
- Conditioning buffer	Phosphate 0.1M pH7 + 0.1 % BSA
- Appearance	Frozen form
- Storage conditions	<-16 °C
- Expiration date	2030.05.18
- Volume per vial	1.6 mL
- Active Moiety Concentration (Mab)	0.25 mg/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 Life Sciences web site: www.cisbio.com