

**Anti HA - d2**

MAb Anti HA-d2 - 5,000 tests

Product reference :

**610HADAA**

**Product characteristics :**

- Lot Number	06A
- Conditioning buffer	Phosphate 0.1M pH7 + 0.1 % BSA
- Appearance	Frozen form
- Storage conditions	-20 °C
- Expiration date	2027.06.19
- Volume per vial	0.4 mL
- Active Moiety Concentration (Mab)	0.25 mg/mL

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (Mab)	0.10 mg	150 000

*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)*