

**MAb MBP-D2**

MAb Anti MBP-d2 - 5,000 tests

Product reference :

**61MBPDAA**

**Product characteristics :**

- Lot Number	06A
- Packaging Unit	5 000 tests/vial
- Conditioning buffer	Phosphate 0.1M pH7 + 0.1% BSA
- Appearance	Frozen form
- Storage conditions	<-16 °C
- Expiration date	2028.05.27
- Volume per vial	0.4 mL
- Active Moiety Concentration (Mab)	250 µg/mL

**Weight indications :**

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

**Recommended quantity per well\* :**

	in 384 well low volume plate (20µL)
Active Moiety (Mab)	20.00 ng
Number of tests/vial	5 000

\*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)**