

PAb Anti Human IgG-d2

PAb Anti Human IgG-d2 - 5,000 tests

Product reference :

61HFCDA

Product characteristics :

- Lot Number _____ 04B
- Packaging Unit _____ 5 000 tests/vial
- Conditioning buffer _____ Phosphate 0.1M pH7 + 0.1% BSA
- Appearance _____ Frozen form
- Storage conditions _____ <-16 °C
- Expiration date _____ 2024.04.27
- Volume per vial _____ 0.4 mL
- Active Moiety Concentration (PAb) _____ 250 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (PAb)	100.00 µg	150 000

Recommended quantity per well* :

	in 384 well low volume plate (20µL)
Active Moiety (PAb)	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