

PAb Anti Human IgG-d2

Product reference :

PAb Anti Human IgG-d2 - 20,000 tests

61HFCDAB**Product characteristics :**

- Lot Number	06A
- Packaging Unit	20 000 tests/vial
- Conditioning buffer	Phosphate 0.1M pH7 + 0.1% BSA
- Appearance	Frozen form
- Storage conditions	<-16 °C
- Expiration date	2027.12.19
- Volume per vial	1.6 mL
- Active Moiety Concentration (PAb)	250 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (PAb)	0.40 mg	150 000

Recommended quantity per well* :

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