



Technical Data Sheet

PAb Anti Human IgG-d2

HTRF PAb Anti Human IgG-d2 - 20K pts

Product reference :

61HFCDAB

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 2034.01.29
- Volume per vial 1.60 mL
- Active Moiety Concentration (PAb) 250.00 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (PAb)	400.00 µg	150 000

Recommended quantity per well* :

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

*Actual amount per well should be optimized according to assay conditions

PLEASE NOTE

Further information as well as a specific application note regarding conjugate-related calculations are available on our Life Sciences web site: www.cisbio.com