

PAb Anti human IgG-XL

Product reference :

PAb Anti Human IgG-XL665 - 5,000 tests

61HFCXLA**Product characteristics :**

- Lot Number	13D
- Packaging Unit	5 000 tests/vial
- Conditioning buffer	Phosphate 0.1M pH7 + 0.1% BSA + stabilizers
- Appearance	Lyophilized form
- Storage conditions	2-8 °C
- Expiration date	2022.07.15
- Reconstitute each vial with	0.25 mL distilled H2O
- Active Moiety Concentration (PAb)	400 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

PLEASE NOTE

After reconstitution the concentration of Active Moiety (PAb) is 400 µg/mL.

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