

Batch information

ABDNP-XL

MAb Anti DNP-XL665 - 5,000 tests

Product reference :

61DNPXLA

Product characteristics :

- Lot Number	16Y
- 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	2023.10.24
- Reconstitute each vial with	0.25 mL distilled H2O
- Active Moiety Concentration (Ab)	400.00 µg/mL

Weight indications :

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

Recommended quantity per well* :

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

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

PLEASE NOTE

The expiration date of batch 16Y has been extended to 2023-10-24 after validation by QC.

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

