

Batch information

ABDNP-XL

MAb Anti DNP-XL665 - 20,000 tests

Product reference :

61DNPXLB

Product characteristics :

- Lot Number	16X
- Packaging Unit	20 000 tests/vial
- Conditioning buffer	Phosphate 0.1M pH7 + 0.1% BSA + stabilizers
- Appearance	Lyophilized form
- Storage conditions	2-8 °C
- Expiration date	2024.03.18
- Reconstitute each vial with	1 mL distilled H2O

Weight indications :

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

Recommended quantity per well* :

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

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

PLEASE NOTE

The expiration date of batch 16X has been extended to 2024-03-18 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

