

Ab HA-XL

MAb Anti HA-XL665 - 20,000 tests

Product reference :

610HAXLB

Product characteristics :

- Lot Number_____ 13A
- 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_____ 2022.12.04
- Volume per vial_____ 1 mL
- Active Moiety Concentration (Ab)_____ 400 µg/mL

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 13A has been extended to 2022-12-04 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