

**Ab HA-XL**

MAb Anti HA-XL665 - 1,000 tests

Product reference :

**610HAXLF**

**Product characteristics :**

- Lot Number\_\_\_\_\_ 13A
- Packaging Unit\_\_\_\_\_ 1 000 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\_\_\_\_\_ 0.25 mL
- Active Moiety Concentration (Ab)\_\_\_\_\_ 80 µg/mL

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (Ab)	20.00 µg	150 000

**Recommended quantity per well\* :**

	in 384 well low volume plate (20µL)
Active Moiety (Ab)	20.00 ng
Number of tests/vial	1 000

\*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](http://www.cisbio.com)**