

## Batch information

**Ab HA-XL**

HTRF MAb Anti HA-XL665 - 1K pts

Product reference :

**610HAXLF**

### Product characteristics :

- Lot Number	13 A
- 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	2024.12.04
- Volume per vial	0.25 mL
- Active Moiety Concentration (Ab)	80.00 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 13 A has been extended to 2024-12-04 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)**

