

## Batch information

**Ab HA-XL**

HTRF MAb Anti HA-XL665 - 20K pts

Product reference :

**610HAXLB**

### Product characteristics :

- Lot Number	15A
- 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	2024.11.08
- Reconstitute each vial with	1 mL distilled H2O
- Active Moiety Concentration (Ab)	400.00 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

**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)**

