

## Batch information

**Ab HA-XL**

MAb Anti HA-XL665 - 5,000 tests

Product reference :

**610HAXLA**

### Product characteristics :

|                                    |  |
|------------------------------------|--|
| - Lot Number                       | 14A  |
| - Packaging Unit                   | 5 000 tests/vial                             |
| - Conditioning buffer              | Phosphate 0.1M pH7 + 0.1 % BSA + stabilizers |
| - Appearance                       | Lyophilized form                             |
| - Storage conditions               | 2-8 °C                                       |
| - Expiration date                  | 2023.11.20                                   |
| - Reconstitute each vial with      | 0.25 mL distilled H2O                        |
| - Active Moiety Concentration (Ab) | 400.00 µg/mL                                 |

### Weight indications :

|                    | Quantity per vial | MW (Da) |
|--------------------|-------------------|---------|
| Active Moiety (Ab) | 100.00 µg         | 150 000 |

### Recommended quantity per well\* :

|                      | in 384 well low volume plate (20µL) |
|----------------------|-------------------------------------|
| Active Moiety (Ab)   | 20.00 ng                            |
| Number of tests/vial | 5 000                               |

\*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)**

