

## Batch information

**SA-XL**

Streptavidin-XL665 - 20,000 tests

Product reference :

**610SAXLB**

### Product characteristics :

|                                    |  |
|------------------------------------|--|
| - Lot Number                       | 166A   |
| - 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                  | 2023.12.16                                   |
| - Reconstitute each vial with      | 1 mL distilled H2O                           |
| - Active Moiety Concentration (SA) | 1.00 mg/mL                                   |

### Weight indications :

|                    | Quantity per vial | MW (Da) |
|--------------------|-------------------|---------|
| Active Moiety (SA) | 1.00 mg           | 60 000  |

### Recommended quantity per well\* :

|                      | in 384 well low volume plate (20µL) |
|----------------------|-------------------------------------|
| Active Moiety (SA)   | 50.00 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)**

