

Batch information

SA-XL

Streptavidin-XL665 - 100,000 tests

Product reference :

610SAXAC

Product characteristics :

| | |
|------------------------------------|--|
| - Lot Number | 167C |
| - Packaging Unit | 100 000 tests/vial |
| - Conditioning buffer | Phosphate 0.1M pH7 + 0.1 % BSA + stabilizers |
| - Appearance | Frozen form |
| - Storage conditions | <-60 °C |
| - Expiration date | 2031.04.20 |
| - Volume per vial | 5.00 mL |
| - Active Moiety Concentration (SA) | 1.00 mg/mL |

Weight indications :

| | Quantity per vial | MW (Da) |
|--------------------|-------------------|---------|
| Active Moiety (SA) | 5.00 mg | 60 000 |

Recommended quantity per well* :

| | in 384 well low volume plate (20µL) |
|----------------------|-------------------------------------|
| Active Moiety (SA) | 50.00 ng |
| Number of tests/vial | 100 000 |

*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

