

SA-XL

Streptavidin-XL665 - 3 mg

Product reference :

610SAXLG**Product characteristics :**

| | |
|------------------------------------|--|
| - Lot Number | 162A |
| - Packaging Unit | 60 000 tests/vial |
| - Conditioning buffer | Phosphate 0.1M pH7 + 0.1 % BSA + stabilizers |
| - Appearance | Lyophilized form |
| - Storage conditions | 2-8 °C |
| - Expiration date | 2023.07.28 |
| - Reconstitute each vial with | 3 mL distilled H2O |
| - Active Moiety Concentration (SA) | 1 mg/mL |

Weight indications :

| | Quantity per vial | MW (Da) |
|--------------------|-------------------|---------|
| Active Moiety (SA) | 3.00 mg | 60 000 |

Recommended quantity per well* :

| | in 384 well low volume plate (20µL) |
|----------------------|-------------------------------------|
| Active Moiety (SA) | 50.00 ng |
| Number of tests/vial | 60 000 |

*Actual amount per well should be optimized according to assay conditions

Further information as well as a specific application note regarding conjugate-related calculations are available on our Life Sciences web site: www.cisbio.com