

SA-XL

Streptavidin-XL665 - 3 mg

Product reference :

610SAXLG

Product characteristics :

- Lot Number_____ 162B
- 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 |

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

