

SA-XL

Streptavidin-XL665 - 100,000 tests - CUSTOM LABELING

Product reference :

610SAXAC

Product characteristics :

- Lot Number_____ 158B
- Conditioning buffer_____ Phosphate 0.1M pH7 + 0.1 % BSA
- Appearance_____ Frozen form
- Storage conditions_____ <-60 °C
- Expiration date_____ 2030.03.07
- Volume per vial_____ 5 mL
- Active Moiety Concentration (SA)_____ 1 mg/mL

Weight indications :

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