

SA-XL

Streptavidin-XL665 - 100,000 tests - CUSTOM LABELING
610SAXAC

Product reference :

Product characteristics :

- Lot Number _____ 160A
- Packaging Unit _____ 100 000 tests/vial
- Conditioning buffer _____ Phosphate 0.1M pH7 + 0.1 % BSA
- Appearance _____ Frozen form
- Storage conditions _____ <-60 °C
- Expiration date _____ 2030.04.01
- 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

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

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