

SA-XL

Streptavidin-XL665 - 5,000 tests

Product reference :

610SAXLA

Product characteristics :

- Lot Number_____ 164A
- Conditioning buffer_____ Phosphate 0.1M pH7 + 0.1 % BSA + stabilizers
- Appearance_____ Lyophilized form
- Storage conditions_____ 2-8 °C
- Expiration date_____ 2023.10.02
- Reconstitute each vial with_____ 0.25 mL distilled H2O
- Active Moiety Concentration (SA)_____ 1 mg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (SA)	0.25 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