

SA-XLent!

Streptavidin-XLent! - 20,000 tests

Product reference :

611SAXLB

Product characteristics :

- Lot Number_____ 26A
- Conditioning buffer_____ Phosphate 0.1M pH7 + 0.1% BSA + stabilizers
- Appearance_____ Lyophilized form
- Storage conditions_____ 2-8 °C
- Expiration date_____ 2023.05.03
- Reconstitute each vial with_____ 1 mL distilled H2O
- Active Moiety Concentration (SA)_____ 1 mg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (SA)	1.00 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

