

SA-XLent !

Streptavidin-XLent! - 5,000 tests
611SAXLA

Product reference :

Product characteristics :

- Lot Number_____ 25A
- Packaging Unit_____ 5 000 tests/vial
- Conditioning buffer_____ Phosphate 0.1M pH7 + 0.1% BSA + stabilizers
- Appearance_____ Lyophilized form
- Storage conditions_____ 2-8 °C
- Expiration date_____ 2022.06.26
- 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)	250.00 µg	60 000

Recommended quantity per well* :

	in 384 well low volume plate (20µL)
Active Moiety (SA)	50.00 ng
Number of tests/vial	5 000

*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

