

SA-XL

Streptavidin-XL665 - 20,000 tests

Product reference :

610SAXLB**Product characteristics :**

- Lot Number	161A
- Packaging Unit	20 000 tests/vial
- Conditioning buffer	Phosphate 0.1M pH7 + 0.1 % BSA + stabilizers
- Appearance	Lyophilized form
- Storage conditions	2-8 °C
- Expiration date	2023.05.14
- 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

Recommended quantity per well* :

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

*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