

SA-XL

Streptavidin-XL665 - 5,000 tests

Product reference :

610SAXLA**Product characteristics :**

- Lot Number	151A
- 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	2022.07.24
- Reconstitute each vial with	250 µL 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