

SA-XL

Streptavidin-XL665 - 3 mg

Product reference :

610SAXLG

Product characteristics :

- Lot Number_____ 155A
- Packaging Unit_____ 60 000 tests/vial
- Conditioning buffer_____ Phosphate 0.1M pH7 + 0.1 % BSA + stabilizers
- Appearance_____ Lyophilized form
- Storage conditions_____ 2-8 °C
- Expiration date_____ 2022.12.11
- Reconstitute each vial with_____ 3 mL distilled H2O
- Active Moiety Concentration (SA)_____ 1 mg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (SA)	3.00 mg	60 000

Recommended quantity per well* :

	in 384 well low volume plate (20µL)
Active Moiety (SA)	50.00 ng
Number of tests/vial	60 000

*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