

**SA-XL**

Streptavidin-XL665 - 3 mg

Product reference :

**610SAXLG****Product characteristics :**

- Lot Number	148A
- 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	2022.02.06
- 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](http://www.cisbio.com)**