

**SA-XL**

Streptavidin-XL665 - 3 mg

Product reference :

**610SAXLG**

**Product characteristics :**

- Lot Number\_\_\_\_\_ 158B
- Conditioning buffer\_\_\_\_\_ Phosphate 0.1M pH7 + 0.1 % BSA + stabilizers
- Appearance\_\_\_\_\_ Lyophilized form
- Storage conditions\_\_\_\_\_ 2-8 °C
- Expiration date\_\_\_\_\_ 2023.02.12
- 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

**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)**