

**SA-XL**

Streptavidin-XL665 - 3 mg

Product reference :

**610SAXLG**

**Product characteristics :**

- Lot Number\_\_\_\_\_ 159A  
- Conditioning buffer\_\_\_\_\_ Phosphate 0.1M pH7 + 0.1 % BSA + stabilizers  
- Appearance\_\_\_\_\_ Lyophilized form  
- Storage conditions\_\_\_\_\_ 2-8 °C  
- Expiration date\_\_\_\_\_ 2023.03.08  
- 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)**