

## Batch information

**SA-XL** Streptavidin-XL665 - 3 mg  
Product reference : **610SAXLG**

### Product characteristics :

- Lot Number	169A
- 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	2024.06.23
- Reconstitute each vial with	3 mL distilled H2O
- Active Moiety Concentration (SA)	1.00 mg/mL

### Weight indications :

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

### Recommended quantity per well\* :

	in Test (384well lo)
Active Moiety (SA)	50.00 ng
Number of tests/vial	60 000

\*Actual amount per well should be optimized according to assay conditions

### PLEASE NOTE

**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)**

