

**SA-XL**

Streptavidin-XL665 - 5,000 tests

Product reference :

**610SAXLA**

**Product characteristics :**

- Lot Number\_\_\_\_\_ 158A
- Packaging Unit\_\_\_\_\_ 5 000 tests/vial
- 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\_\_\_\_\_ 0.25 mL distilled H2O
- Active Moiety Concentration (SA)\_\_\_\_\_ 1 mg/mL

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (SA)	250.00 µg	60 000

**Recommended quantity per well\* :**

	in 384 well low volume plate (20µL)
Active Moiety (SA)	50.00 ng
Number of tests/vial	5 000

\*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)**