

**SA-XL**

Streptavidin-XL665 - 100,000 tests - CUSTOM LABELING  
**610SAXAC**

Product reference :

**Product characteristics :**

- Lot Number \_\_\_\_\_ 137A
- Packaging Unit \_\_\_\_\_ 100 000 tests/vial
- Conditioning buffer \_\_\_\_\_ Phosphate 0.1M pH7 + 0.1 % BSA
- Appearance \_\_\_\_\_ Frozen form
- Storage conditions \_\_\_\_\_ 2-8 °C
- Expiration date \_\_\_\_\_ 2026.10.22
- Volume per vial \_\_\_\_\_ 6.25 mL
- Active Moiety Concentration (SA) \_\_\_\_\_ 0.8 mg/mL

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (SA)	5.00 mg	60 000

**Recommended quantity per well\* :**

	in 384 well low volume plate (20µL)
Active Moiety (SA)	50.00 ng
Number of tests/vial	100 000

\*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)**

