

**SA-XL**

Streptavidin-XL665 - 100,000 tests

Product reference :

**610SAXAC**

**Product characteristics :**

- Lot Number\_\_\_\_\_ 134A
- Packaging Unit\_\_\_\_\_ 100 000 tests/vial
- Conditioning buffer\_\_\_\_\_ Phosphate 0.1M pH7 + 0.1 % BSA + stabilizers
- Appearance\_\_\_\_\_ Frozen form
- Storage conditions\_\_\_\_\_ <-60 °C
- Expiration date\_\_\_\_\_ 2025.09.14
- Volume per vial\_\_\_\_\_ 5 mL
- Active Moiety Concentration (SA)\_\_\_\_\_ 1 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 HTRF web site: [www.htrf.com/products](http://www.htrf.com/products)**