

Batch information

SA-XL

CUSTOM LABELING

Product reference :

610SAXAC

Product characteristics :

- Lot Number	170A
- Packaging Unit	100 000 tests/vial
- Conditioning buffer	Phosphate 0.1M pH7 + 0.1 % BSA
- Appearance	Frozen form
- Storage conditions	<-60° °C
- Expiration date	2031.07.07
- Volume per vial	5.00 mL
- Active Moiety Concentration (SA)	1.00 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

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

