

Batch information

SA-XL

CUSTOM LABELING

Product reference :

610SAXAC

Product characteristics :

- Lot Number	172A
- Conditioning buffer	Phosphate 0.1M pH7 + 0.1 % BSA
- Appearance	Frozen form
- Storage conditions	<-60° °C
- Expiration date	2031.11.08
- 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

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

