

Batch information

STD Insulin kit - Standard

Product reference : **62INSCDA**

Product characteristics :

- Lot Number	16A
- Conditioning buffer	Phosphate 50 mM pH7.0 + BSA 0.2% + stabilizers
- Appearance	Lyophilized form
- Storage conditions	2-8 °C
- Expiration date	2024.06.02

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (ins)	225.00 ng	5 807

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

