

STD Insulin kit calibrator
Product reference : **62INSCDA**

Product characteristics :

- Lot Number_____ 14A
- Conditioning buffer_____ Phosphate 50 mM pH7.0 +
BSA 0.2% + stabilizers
- Appearance_____ Lyophilized form
- Storage conditions_____ 2-8 °C
- Expiration date_____ 2022.06.08

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (ins)	275.00 ng	5 807

PLEASE NOTE

The expiration date of batch 14A has been extended to 2022-06-08 after validation by QC

Further information as well as a specific application note regarding conjugate-related calculations are available on our Life Sciences web site: www.cisbio.com

