

**STD**

Insulin kit calibrator

Product reference :

**62INSCDA**

**Product characteristics :**

- Lot Number\_\_\_\_\_ 15A  
- Conditioning buffer\_\_\_\_\_ Phosphate 50 mM pH7.0 +  
BSA 0.2% + stabilizers  
- Appearance\_\_\_\_\_ Lyophilized form  
- Storage conditions\_\_\_\_\_ 2-8 °C  
- Expiration date\_\_\_\_\_ 2023.02.12

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (ins)	200.00 ng	5 807

*Further information as well as a specific application note regarding conjugate-related calculations are available on our Life Sciences web site: [www.cisbio.com](http://www.cisbio.com)*