

**Rec buffer 1**

Reconstitution buffer 1 - 16 ml

Product reference :

**62RB1RDD**

**Product characteristics :**

- Lot Number \_\_\_\_\_ 12A  
- Conditioning buffer \_\_\_\_\_ -  
- Appearance \_\_\_\_\_ Liquid form  
- Storage conditions \_\_\_\_\_ 2-8 °C  
- Expiration date \_\_\_\_\_ 2024.11.11

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (Buf)	16.00 mL	N/A

*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)*