

**Binding Domain diluent buffer**      Binding Domain diluent buffer - 200 ml  
Product reference :                      **62DLBDDF**

**Product characteristics :**

- Lot Number \_\_\_\_\_ 07A  
- Conditioning buffer \_\_\_\_\_ -  
- Appearance \_\_\_\_\_ Liquid form  
- Storage conditions \_\_\_\_\_ -20 °C  
- Expiration date \_\_\_\_\_ 2023.01.30  
- Volume per vial \_\_\_\_\_ 200 mL

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (Buf)	200.00 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)*