

**Diluent 3**

Product reference :

Diluent 3 - 20 ml

**62DL3DDD**

**Product characteristics :**

- Lot Number \_\_\_\_\_ 10A  
- Conditioning buffer \_\_\_\_\_ -  
- Appearance \_\_\_\_\_ Liquid form  
- Storage conditions \_\_\_\_\_ 2-8 °C  
- Expiration date \_\_\_\_\_ 2023.06.04

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (Buf)	20.00 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)*