

**Diluent 2**

Product reference :

Diluent 2 - 20 mL

**62DL2DDD**

**Product characteristics :**

- Lot Number\_\_\_\_\_ 09A  
- Conditioning buffer\_\_\_\_\_ -  
- Appearance\_\_\_\_\_ Liquid form  
- Storage conditions\_\_\_\_\_ 2-8 °C  
- Expiration date\_\_\_\_\_ 2024.02.11

**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)*