

**Diluent 5**

Product reference :

Diluent 5 (5X) - 2 ml

**62DL5DDA**

**Product characteristics :**

- Lot Number \_\_\_\_\_ 16C  
- Conditioning buffer \_\_\_\_\_ -  
- Appearance \_\_\_\_\_ Frozen form  
- Storage conditions \_\_\_\_\_ <-16 °C  
- Expiration date \_\_\_\_\_ 2023.06.19

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (Buf)	2.00 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)*

