

**Diluent 9**

Product reference :

Diluent 9 (5X) - 2 ml

**62DL9DDA**

**Product characteristics :**

- Lot Number\_\_\_\_\_ 01C  
- Conditioning buffer\_\_\_\_\_ -  
- Appearance\_\_\_\_\_ Frozen form  
- Storage conditions\_\_\_\_\_ <-16 °C  
- Expiration date\_\_\_\_\_ 2024.06.30

**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)*