

Diluent 5

Product reference :

Diluent 5 (5X) - 2 ml

62DL5DDA

Product characteristics :

- Lot Number _____ 15C
- Conditioning buffer _____ -
- Appearance _____ Frozen form
- Storage conditions _____ <-16 °C
- Expiration date _____ 2023.01.22

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

