

Diluent 5

Product reference :

Diluent 5 (5X) - 100 ml

62DL5DDG

Product characteristics :

- Lot Number _____ 20A
- Conditioning buffer _____ -
- Appearance _____ Liquid form
- Storage conditions _____ 2-8 °C
- Expiration date _____ 2025.03.12

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Buf)	100.00 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

