

Diluent 9

Product reference :

Diluent 9 (5X) - 100 ml

62DL9DDC

Product characteristics :

- Lot Number_____ 02A
- Conditioning buffer_____ -
- Appearance_____ Liquid form
- Storage conditions_____ 2-8 °C
- Expiration date_____ 2024.05.26

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