

Diluent 5-10ml
Product reference :

Diluent 5 (5X) - 10 ml
62DL5DDC

Product characteristics :

- Lot Number _____ 18A
- Conditioning buffer _____ -
- Appearance _____ Liquid form
- Storage conditions _____ 2-8 °C
- Expiration date _____ 2024.11.23

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Buf)	10.00 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