

Rec buffer 3

Reconstitution buffer 3 - 200 mL

Product reference :

62RB3RDF

Product characteristics :

- Lot Number _____ 23A
- Conditioning buffer _____ -
- Appearance _____ Liquid form
- Storage conditions _____ 2-8 °C
- Expiration date _____ 2023.12.06

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Buf)	200.00 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

