

Elution Buffer

Product reference :

Elution buffer 1 - 20 ml

62EB1FDD

Product characteristics :

- Lot Number_____ 13A
- Conditioning buffer_____ -
- Appearance_____ Liquid form
- Storage conditions_____ 2-8 °C
- Expiration date_____ 2022.11.14
- Volume per vial_____ 20 mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Buf)	20.00 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