

Batch information

Stimulation buffer #4 Stimulation buffer 4 (5X) - 200 mL
Product reference : **64SB4FDF**

Product characteristics :

- Lot Number 15A
- Conditioning buffer -
- Appearance Liquid form
- Storage conditions 2-8 °C
- Expiration date 2026.03.19

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Tp)	200.00 mL	N/A

PLEASE NOTE

Further information as well as a specific application note regarding conjugate-related calculations are available on our Life Sciences web site: www.cisbio.com

