

Stimulation buf
Product reference :

Stimulation Buffer 2 (5X) - 8 ml
62IP1FDC

Product characteristics :

- Lot Number _____ 32A
- Conditioning buffer _____ -
- Appearance _____ Liquid form
- Storage conditions _____ 2-8 °C
- Expiration date _____ 2023.07.10

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Tp)	8.00 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

