

Batch information

Diluent 5 HTRF Diluent 5 (5X) - 100 mL
Product reference : **62DL5DDG**

Product characteristics :

- Lot Number 21A
- Conditioning buffer -
- Appearance Liquid form
- Storage conditions 2-8 °C
- Expiration date 2028.10.04

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Buf)	100.00 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

