

Batch information

Diluent 3 HTRF Diluent 3 - 20 mL
Product reference : **62DL3DDD**

Product characteristics :

- Lot Number 12A
- Conditioning buffer -
- Appearance Liquid form
- Storage conditions 2-8 °C
- Expiration date 2026.10.28

Weight indications :

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

