

Batch information

Diluent 5-10ml

HTRF Diluent 5 (5X) - 10 mL

Product reference :

62DL5DDC

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)	10.00 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

