

Batch information

Binding Domain Detection buffer 1 HTRF Bind. Domain - Detect. Buf.1-130 mL

Product reference : **62DB1FDG**

Product characteristics :

- Lot Number	11A
- Conditioning buffer	-
- Appearance	Liquid form
- Storage conditions	2-8 °C
- Expiration date	2024.05.26
- Volume per vial	130.00 mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Buf)	130.00 mL	1

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

