

## Batch information

**Binding Domain diluent buffer** Binding Domain - Diluent buffer- 200 mL

Product reference : **62DLBDDF**

### Product characteristics :

- Lot Number	09A
- Conditioning buffer	-
- Appearance	Liquid form
- Storage conditions	2-8 °C
- Expiration date	2024.07.01
- Volume per vial	200.00 mL

### Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Buf)	200.00 mL	N/A

**Further information as well as a specific application note regarding conjugate-related calculations are available on our Life Sciences web site: [www.cisbio.com](http://www.cisbio.com)**

