

## Batch information

**Diluent 5** HTRF Diluent 5 (5X) - 2 mL  
Product reference : **62DL5DDA**

### Product characteristics :

- Lot Number 19G  
- Conditioning buffer -  
- Appearance Frozen form  
- Storage conditions <-16 °C  
- Expiration date 2025.01.22

### Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Buf)	2.00 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](http://www.cisbio.com)*

