

## Batch information

**Detection buffer** HTRF PPI-Terbium Detect. Buf. - 220 mL  
Product reference : **61DB10RDF**

### Product characteristics :

- Lot Number 08A  
- Conditioning buffer -  
- Appearance Liquid form  
- Storage conditions 2-8 °C  
- Expiration date 2027.03.27

### Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Buf)	220.00 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)*

