

**Detection buffer**

PPI - Europium detection buffer - 220 mL

Product reference :

**61DB9RDF**

**Product characteristics :**

- Lot Number \_\_\_\_\_ 03A  
- Conditioning buffer \_\_\_\_\_ -  
- Appearance \_\_\_\_\_ Liquid form  
- Storage conditions \_\_\_\_\_ 2-8 °C  
- Expiration date \_\_\_\_\_ 2024.04.08

**Weight indications :**

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