

Batch information

Detection buffer HTRF PPI-Europium Detect. Buf. - 220 mL
Product reference : **61DB9RDF**

Product characteristics :

- Lot Number 07A
- Conditioning buffer -
- Appearance Liquid form
- Storage conditions 2-8 °C
- Expiration date 2027.02.09

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

