

## Batch information

**EPiGeneous detection buffer One** HTRF Epig. Detect. Buf. 1 - 24 mL

Product reference : **62SAHRDD**

### Product characteristics :

- Lot Number 15A  
- Conditioning buffer -  
- Appearance Frozen form  
- Storage conditions ==-20 °C  
- Expiration date 2024.10.09  
- Volume per vial 24.00 mL

### Weight indications :

|                     | Quantity per vial | MW (Da) |
|---------------------|-------------------|---------|
| Active Moiety (Buf) | 24.00 mL          | N/A     |

### PLEASE NOTE

The expiration date of batch 15A has been extended to 2024-10-09 after validation by QC.

***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)***

