

## Batch information

### Stabilization buffer 1

HTRF Stabilization Buf. 1 - 200 mL

Product reference :

**62SB1FDF**

### Product characteristics :

- Lot Number	03A
- Conditioning buffer	-
- Appearance	Frozen form
- Storage conditions	2-8 °C
- Expiration date	2027.06.16

### Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (buf)	200.00 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)*

