

**Stimulation Buf**  
Product reference :Stimulation Buffer 1 (5X) - 8 ml  
**64SB1FDC****Product characteristics :**

- Lot Number \_\_\_\_\_ 08A  
- Conditioning buffer \_\_\_\_\_ -  
- Appearance \_\_\_\_\_ Liquid form  
- Storage conditions \_\_\_\_\_ 2-8 °C  
- Expiration date \_\_\_\_\_ 2023.11.20

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (Tp)	8.00 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)*