

Binding Domain Detection buffer #1 Binding Domain Detection buffer #1 - 130 ml
Product reference : **62DB1FDG**

Product characteristics :

- Lot Number _____ 10A
- Conditioning buffer _____ -
- Appearance _____ Frozen form
- Storage conditions _____ -20 °C
- Expiration date _____ 2023.07.10
- Volume per vial _____ 130 mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Buf)	130.00 mL	1

Further information as well as a specific application note regarding conjugate-related calculations are available on our Life Sciences web site: www.cisbio.com