

EPIgeneous detection buffer One EPIgeneous detection buffer one - 24 ml
Product reference : **62SAHRDD**

Product characteristics :

- Lot Number_____ 15A
- Conditioning buffer_____ -
- Appearance_____ Frozen form
- Storage conditions_____ =-20 °C
- Expiration date_____ 2022.10.09
- Volume per vial_____ 24 mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Buf)	24.00 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