

EPIgeneous detection buffer Two EPIgeneous detection buffer two - 130 ml
Product reference : **62SAHFDD**

Product characteristics :

- Lot Number_____ 12A
- Conditioning buffer_____ -
- Appearance_____ Frozen form
- Storage conditions_____ < -20 °C
- Expiration date_____ 2023.12.12
- Volume per vial_____ 130 mL

Weight indications :

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