

Batch information

EPi geneous detection buffer Two HTRF Epig. Detect. Buf. 2 - 130 mL

Product reference : **62SAHFDD**

Product characteristics :

- Lot Number 13A
- Conditioning buffer -
- Appearance Frozen form
- Storage conditions < -20 °C
- Expiration date 2027.04.18
- Volume per vial 130.00 mL

Weight indications :

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

