

Batch information

Control lysate HTRF S1/ACE2 Bind. kit - Ctrl
Product reference : **64BDS1TDA**

Product characteristics :

- Lot Number 01A
- Conditioning buffer Tp PO4 50mM + KF 0.8M + 0.2% BSA pH7
- Appearance Frozen form
- Storage conditions -60 °C
- Expiration date 2023.10.20

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (0)	50.00 µl	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

