

Batch information

Control lysate HTRF SHP1 total kit - Ctrl lysate
Product reference : **64NH1TDA**

Product characteristics :

- Lot Number 01A
- Conditioning buffer lysis buffer
- Appearance Frozen form
- Storage conditions <-60 °C
- Expiration date 2023.07.03
- Volume per vial 150.00 µL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (mAb)	150.00 µl	N/A

PLEASE NOTE

The expiration date of batch 01A has been extended to 2023-07-03 after validation by QC.

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

