



Technical Data Sheet

Control lysate

HTRF (h) CHK2 p-T68 Kit - Ctrl Lysate

Product reference :

64CHK2T8TDA

Product characteristics :

- Lot Number 02AR
- Conditioning buffer Lysis Buffer
- Appearance Frozen form
- Storage conditions -60 °C
- Expiration date 2026.11.28
- Volume per vial 150.00 µL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (CL)	150.00 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