



Technical Data Sheet

Tau Total control lysate

HTRF Tau Total kit - Ctrl lysate

Product reference :

64NTAUTDA

Product characteristics :

- Lot Number	04AR
- Conditioning buffer	Cellular kinase lysis buffer
- Appearance	Frozen form
- Storage conditions	-60 °C
- Expiration date	2027.01.02

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