

Control lysate

ERK total kit control lysate

Product reference :

64NRKTDA

Product characteristics :

- Lot Number _____ 10A
- Conditioning buffer _____ Cellular kinase lysis buffer
- Appearance _____ Frozen form
- Storage conditions _____ -60 °C
- Expiration date _____ 2022.04.02
- Volume per vial _____ 150 µL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Ab)	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

