

Batch information

Control lysate

STING total kit - control lysate

Product reference :

64NTGTDA

Product characteristics :

- Lot Number	05A
- Conditioning buffer	Phospho total protein lysis buffer
- Appearance	Frozen form
- Storage conditions	-60 °C
- Expiration date	2024.06.08
- Volume per vial	150.00 µL
- Active Moiety Concentration (Ab)	1.00 X

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

