

Control lysate

SLP-76 total control lysate

Product reference :

63ADK077TDA

Product characteristics :

- Lot Number	03A
- Conditioning buffer	Lysis Buffer #1
- Appearance	Frozen form
- Storage conditions	2-8 °C
- Expiration date	2023.01.09
- Volume per vial	150 µL

Weight indications :

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

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