

Control lysate

SLP-76 phospho S376 control lysate

Product reference :

63ADK076TDA

Product characteristics :

- Lot Number _____ 04A
- Conditioning buffer _____ Lysis Buffer # 1
- Appearance _____ Frozen form
- Storage conditions _____ -60 °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