

Control lysate

FoxO1 phospho-S256 kit control lysate

Product reference :

64FOXTDA

Product characteristics :

- Lot Number _____ 01A
- Conditioning buffer _____ Lysis Buffer
- Appearance _____ Frozen form
- Storage conditions _____ -60 °C
- Expiration date _____ 2023.01.18
- Volume per vial _____ 150 µL

Weight indications :

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

PLEASE NOTE

The expiration date of batch 01A has been extended to 2023-01-18 after validation by QC

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

