

Control lysate

Beta Catenin phospho-T41/ S33/37 control lysate

Product reference :

64CATTDA

Product characteristics :

- Lot Number_____ 02A
- Conditioning buffer_____ Lysis Buffer
- Appearance_____ Frozen form
- Storage conditions_____ -60 °C
- Expiration date_____ 2022.11.07
- 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 02A has been extended to 2022-11-07 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

