

Control lysate

Product reference :

IKKB phospho-S177/181 kit control lysate
64KKBTDA

Product characteristics :

- Lot Number _____ 06A
- Conditioning buffer _____ Phospho total protein lysis buffer
- Appearance _____ Frozen form
- Storage conditions _____ -60 °C
- Expiration date _____ 2022.02.21
- Volume per vial _____ 150 µL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (IKK)	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