

AKT phospho T308 kit control lysate AKT phospho-T308 kit control lysate
Product reference : **62AKTTDA**

Product characteristics :

- Lot Number _____ 06A
- Conditioning buffer _____ Phospho total protein lysis buffer
- Appearance _____ Frozen form
- Storage conditions _____ -60 °C
- Expiration date _____ 2022.02.18
- Volume per vial _____ 150 µL

Weight indications :

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

