

AKT phospho S473 control lysate AKT phospho-S473 kit control lysate
Product reference : **62AKSTDA**

Product characteristics :

- Lot Number_____ 14A
- Conditioning buffer_____ Phospho total protein lysis buffer
- Appearance_____ Frozen form
- Storage conditions_____ -60 °C
- Expiration date_____ 2023.01.09
- Volume per vial_____ 150 µL
- Active Moiety Concentration (Ab)_____ 1 X

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

