

Batch information

AKT phospho S473 control lysate AKT phospho-S473 kit - control lysate

Product reference : **62AKSTDA**

Product characteristics :

- Lot Number	16A
- Conditioning buffer	Phospho total protein lysis buffer
- Appearance	Frozen form
- Storage conditions	-60 °C
- Expiration date	2024.02.16
- Volume per vial	150.00 µL
- Active Moiety Concentration (Ab)	1.00 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

