

## Batch information

**AKT phospho S473 control lysate** HTRF AKT p-S473 kit - Ctrl lysate

Product reference : **62AKSTDA**

### Product characteristics :

- Lot Number	17A
- Conditioning buffer	Phospho total protein lysis buffer
- Appearance	Frozen form
- Storage conditions	-60 °C
- Expiration date	2024.09.30
- 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](http://www.cisbio.com)*

