

## Batch information

### Control lysate

FAK phospho-Y397 kit - control lysate

Product reference :

**64FAKTDA**

### Product characteristics :

- Lot Number	03A
- Conditioning buffer	lysis buffer
- Appearance	Frozen form
- Storage conditions	-60 °C
- Expiration date	2024.10.27
- Volume per vial	150.00 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](http://www.cisbio.com)*

