

## Batch information

**Control lysate** FGFR3 p-Y647/648 kit - control lysate  
Product reference : **64FGFR3Y6TDA**

### Product characteristics :

- Lot Number	01A
- Conditioning buffer	Lysis buffer
- Appearance	Frozen form
- Storage conditions	<-60 °C
- Expiration date	2023.09.22

### 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)*

