

## Batch information

**Control lysate**

FGFR3 p-Y647/648 kit - control lysate

Product reference :

**64FGFR3Y6TDA****Product characteristics :**

- Lot Number	03A
- Conditioning buffer	Lysis buffer
- Appearance	Frozen form
- Storage conditions	<-60 °C
- Expiration date	2024.09.28

**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)***

