

Batch information

control lysate

FGFR1 p-Y653/654 kit - control lysate

Product reference :

64FGFR1Y6TDA

Product characteristics :

- Lot Number	02A
- Conditioning buffer	Lysis buffer
- Appearance	Frozen form
- Storage conditions	<-60 °C
- Expiration date	2024.03.31

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (CL)	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

