

Batch information

Control lysate

FGFR4 p-Y642 kit - control lysate

Product reference :

64FGFR4Y6TDA

Product characteristics :

- Lot Number	01A
- Conditioning buffer	Lysis buffer
- Appearance	Frozen form
- Storage conditions	<-60 °C
- Expiration date	2023.10.01

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

