

Batch information

lysis buf5

Lysis & Detection Buffer 5 - 13 ml

Product reference :

62CL5FDD

Product characteristics :

- Lot Number	06A
- Conditioning buffer	-
- Appearance	Liquid form
- Storage conditions	2-8 °C
- Expiration date	2025.10.22

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Buf)	13.00 mL	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

