

conj/lysis buf1
Product reference :

Lysis & Detection Buffer 1 - 13 ml
62CL1FDD

Product characteristics :

- Lot Number _____ 13A
- Conditioning buffer _____ -
- Appearance _____ Liquid form
- Storage conditions _____ 2-8 °C
- Expiration date _____ 2023.07.04

Weight indications :

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

Further information as well as a specific application note regarding conjugate-related calculations are available on our Life Sciences web site: www.cisbio.com