

Batch information

STD IP-One Gq kit - Standard

Product reference : **62IP1CDA**

Product characteristics :

- Lot Number	13A
- Conditioning buffer	Phosphate 10 mM pH7.0 + BSA 0.2%
- Appearance	Lyophilized form
- Storage conditions	2-8 °C
- Expiration date	2022.07.01

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (IP1)	0.11 µmoles	N/A

PLEASE NOTE

Reconstitute with 0.500 mL H₂O.

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

