

STD

Pro-collagen Type 1 (human) kit- Standard
63ADK014CDA

Product reference :

Product characteristics :

- Lot Number_____ 13A
- Conditioning buffer_____ Phosphate 50 mM pH7.0 +
BSA 0.2% + stabilizers
- Appearance_____ Frozen form
- Storage conditions_____ -80 °C
- Expiration date_____ 2022.01.15

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Ab)	0.50 µg	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