

Batch information

STD Estradiol kit - Standard

Product reference : **62ESTCDA**

Product characteristics :

- Lot Number	05A
- Conditioning buffer	Phosphate 50 mM pH7.0 + BSA 0.2% + stabilizers
- Appearance	Lyophilized form
- Storage conditions	2-8 °C
- Expiration date	2024.06.23

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Ab)	50.00 ng	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

