

Batch information

STD GST check kit - Standard

Product reference : **62GSTCDA**

Product characteristics :

- Lot Number	09A
- Conditioning buffer	Phosphate 50 mM pH7.0 + BSA 0.2% + stabilizers
- Appearance	Lyophilized form
- Storage conditions	2-8 °C
- Expiration date	2023.11.30

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (GST)	0.20 nmoles	N/A

PLEASE NOTE

The expiration date of batch 09A has been extended to 2023-11-30 after validation by QC.

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

