

Product Number:	102335
Lot Number:	125608
Expiration:	13-Sep-2023
Storage:	-20°C

Calibrator Concentrate Target Value (ng/mL)	20
--	----

Laboratory Analysis

Tested with Paired Kit Lot:	n/a
--------------------------------	-----

Review/Approval Sonia Khillan QC Manager Sonia Khillan / 18JAN2023
Name Title Signature/Date