

Product Number:	102335
Lot Number:	108915
Expiration:	30-Mar-2023
Storage:	-20°C

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

Laboratory Analysis

Tested with Paired Kit Lot:	502862
--------------------------------	--------

Review/Approval Sonia Khillan QC Manager Sonia Khillan / 08NOV2022
Name Title Signature/Date