

Product Number:	102335
Lot Number:	312406
Control Number:	312406A
Expiration:	4-May-2025
Storage:	-20°C

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

Laboratory Analysis

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

Review/Approval

ROBERT SOMOGYI	QC MANAGER		03 AUG 2023
Name	Title	Signature/Date	