

Product Number:	103705
Lot Number:	215301
Expiration:	2-Jun-2024
Storage:	-20°C

Laboratory Analysis

Concentration (pg/mL)	974
------------------------------	-----

Review/Approval Sonia Khillan QC Manager Sonia Khillan 07DEC2022
Name Title Signature/Date