

Product Number:	102256
Lot Number:	214303
Expiration:	23-May-2024
Storage:	- 80°C
Kit Lot Number Match:	N/A

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	214303	4.34	3.16 - 4.75

* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

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