

Product Number:	102257
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 2	214303	156	126 - 190

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

Review/Approval

*Sonia Khillan*  
Name Title

*Sanjiv Kalia* 07DEC2022  
Signature/Date