

Product Number:	103650
Lot Number:	301007
Expiration:	10-Apr-2023
Storage:	2-8°C



Data below represents results generated on the Simoa® HD-1 Analyzer.

Laboratory Analysis

Characteristic	Result
HD-1 Analyzer	PASS

SBG Concentrate Concentration (nM)	160
------------------------------------	-----

Review/Approval Sonia Khillan QC Manager Sonia Khillan 102 FEB 2023
 Name Title Signature/Date