

Product Number:	103650
Lot Number:	227803
Expiration:	5-Jan-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
Name

QC Manager
Title

Sonia Khillan / 01DEC2022
Signature/Date