

Product Number:	103718
Lot Number:	503749
Expiration:	5-Jan-2023



Component	Item Number	Lot Number
SBG Concentrate	103650	227803

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

Means of Analysis	Result
HD-1/ HDX	PASS

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

Review/Approval

Sonia Khillan
Name

QC manager
Title

Sonia Rolla / 05 DEC 2022
Signature/Date