

Product Number:	103718
Lot Number:	503800
Expiration:	10-Apr-2023



Component	Item Number	Lot Number
SBG Concentrate	103650	301007

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

Sania Khilkin
Name

QC Manager
Title

Sania Khilkin / 03 FEB 2023
Signature/Date