

Product Number:	103718
Lot Number:	503826
Expiration:	14-May-2023



Component	Item Number	Lot Number
SBG Concentrate	103650	304504

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 Kella 16 MAR 2023
Signature/Date