

Product Number:	103718
Lot Number:	503407
Expiration:	22-May-2022



Component	Item Number	Lot Number
SBG Concentrate	103650	205307

SBG Concentrate Concentration (nM)	138
------------------------------------	-----

Means of Analysis	Result
HD-1/ HDX	PASS

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

Review/Approval

Matthew Smart QC Director
~~Shivani Goel~~ ~~QC Manager~~
 Name W16Mar2022 Title W16Mar2022 Signature/Date MSA 16Mar2022