

Product Number:	103718
Lot Number:	503671
Expiration:	26-Oct-2022



Component	Item Number	Lot Number
SBG Concentrate	103650	220720

SBG Concentrate Concentration (nM)	135
------------------------------------	-----

Means of Analysis	Result
HD-1/ HDX	PASS

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

Review/Approval

Matt Smart
Name

QC
Title

M. A. 15 Aug 2022
Signature/Date