

Product Number:	103718
Lot Number:	503436
Expiration:	15-Jun-2022



Component	Item Number	Lot Number
SBG Concentrate	103650	207404

SBG Concentrate Concentration (nM)	129.8
------------------------------------	-------

Means of Analysis	Result
HD-1/ HDX	PASS

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

Review/Approval

Richa Shrestha
Name

QC Supervisor
Title

[Signature] 07Apr2022
Signature/Date