

Product Number:	103718
Lot Number:	502791
Expiration:	8-May-2021

Component	Item Number	Lot Number
SBG Concentrate	103650	103918

SBG Concentrate Concentration (nM)	139
------------------------------------	-----

Means of Analysis	Result
HD-1/ HDX	PASS

Review/Approval

Junpil Park	QC Supervisor		11Mar2021
Name	Title	Signature/Date	