

Product Number:	103718
Lot Number:	502686
Expiration:	7-Mar-2021

Component	Item Number	Lot Number
SBG Concentrate	103650	034212

SBG Concentrate Concentration (nM)	143
------------------------------------	-----

Means of Analysis	Result
HD-1/ HDX	PASS

Review/Approval

Junpil Park
Name

QC Supervisor
Title



08 Jan 2021

Signature/Date