

Product Number:	103718
Lot Number:	502567
Expiration:	10-Dec-2020

Component	Item Number	Lot Number
SBG Concentrate	103650	025415

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

Means of Analysis	Result
HD-1/ HDX	PASS

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 08 Oct 2020

Name

Title

Signature/Date