

Product Number:	103718
Lot Number:	503064
Expiration:	28-Sep-2021

Component	Item Number	Lot Number
SBG Concentrate	103650	117921

SBG Concentrate Concentration (nM)	152
------------------------------------	-----

Means of Analysis	Result
HD-1/ HDX	PASS

Review/Approval

Shivani Goel

QC Supervisor

Shivani Goel 20 Jul 2021

Name

Title

Signature/Date