

Product Number:	103718
Lot Number:	503296
Expiration:	6-Mar-2022

Component	Item Number	Lot Number
SBG Concentrate	103650	134014

SBG Concentrate Concentration (nM)	128
------------------------------------	-----

Means of Analysis	Result
HD-1/ HDX	PASS

Review/Approval

Shivani Goel  
Name

QC Supervisor  
Title

Shivani Goel 13Jan2022  
Signature/Date