

Product Number:	103718
Lot Number:	503180
Expiration:	27-Dec-2021

Component	Item Number	Lot Number
SBG Concentrate	103650	127010

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

Means of Analysis	Result
HD-1/ HDX	PASS

Review/Approval

Shuvam Goel QC Shuvam Goel ^{SG 13 Oct 2021} _{12 13 Oct 2021}
 Name Title Signature/Date