

Product Number:	103718
Lot Number:	502823
Expiration:	8-Jun-2021

Component	Item Number	Lot Number
SBG Concentrate	103650	106718

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

Means of Analysis	Result
HD-1/ HDX	PASS

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 24Mar2021
Signature/Date