

Product Number:	103718
Lot Number:	502431
Expiration:	4-Sep-2020

Component	Item Number	Lot Number
SBG Concentrate	103650	015605

SBG Concentrate Concentration (nM)	112
------------------------------------	-----

Means of Analysis	Result
HD-1/ HDX	PASS

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 22 Jun 2020
Signature/Date