

Product Number:	103718
Lot Number:	502934
Expiration:	12-Jul-2021

Component	Item Number	Lot Number
SBG Concentrate	103650	110217

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

Means of Analysis	Result
HD-1/ HDX	PASS

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 19 May 2021
Signature/Date