

Product Number:	103718
Lot Number:	502309
Expiration:	26-Jun-2020

Component	Item Number	Lot Number
SBG Concentrate	103650	008606


SBG Concentrate Concentration (nM)	127
------------------------------------	-----

Means of Analysis	Result
HD-1/ HDX	PASS

Review/Approval

Paulo Carvalho  
Name

VP, Manufacturing Operations  
Title

  
Signature/Date

30 Mar 2020  
Date