

Product Number:	101735
Lot Number:	502812
Expiration:	19-Jan-2022

Component	Item Number	Lot Number
SBG Dilution Buffer	100376	104729
SBG Concentrate	103397	101913

Laboratory Analysis

Means of Analysis	Characteristic	Result
HD-1 Analyzer	Functional Test	PASS
SBG Concentrate Concentration (nM)		50

Review/Approval

Junpil Park

Name

QC Supervisor

Title

[Signature] 18 Mar 2021

Signature/Date