

Product Number:	101735
Lot Number:	502834
Expiration:	16-Feb-2022

Component	Item Number	Lot Number
SBG Dilution Buffer	100376	104729
SBG Concentrate	103397	106103

Laboratory Analysis

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

Review/Approval

Jumpil Park
QC Supervisor
[Signature] 26 Mar 2021

Name
Title
Signature/Date