

Product Number:	101735
Lot Number:	502763
Expiration:	8-Jan-2022

Component	Item Number	Lot Number
SBG Dilution Buffer	100376	100805
SBG Concentrate	103397	101912

Laboratory Analysis

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

Review/Approval

Junpil Park

QC Supervisor



22 Feb 2021

Name

Title

Signature/Date