

Product Number:	101735
Lot Number:	502618
Expiration:	27-Aug-2021

Component	Item Number	Lot Number
SBG Dilution Buffer	100376	026623
SBG Concentrate	103397	024007

**Laboratory Analysis**

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

Review/Approval

Shivani Goel  
Name Title

QC Manager

*Shivani Goel* 19 Nov 2020  
Signature/Date