

Product Number:	101735
Lot Number:	502811
Expiration:	15-Feb-2022

Component	Item Number	Lot Number
SBG Dilution Buffer	100376	104729
SBG Concentrate	103397	104620

**Laboratory Analysis**

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

Review/Approval

*Shivani Goel*

Name

*QC Manager*

Title

*Shivani Goel 17 Mar 2021*

Signature/Date