

Product Number:	101735
Lot Number:	502732
Expiration:	16-Oct-2021

Component	Item Number	Lot Number
SBG Dilution Buffer	100376	029015
SBG Concentrate	103397	101116

**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* 08Feb2021  
Signature/Date