

Product Number:	101735
Lot Number:	503209
Expiration:	19-Jul-2022

Component	Item Number	Lot Number
SBG Dilution Buffer	100376	128906
SBG Concentrate	103397	120035

**Laboratory Analysis**

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

Review/Approval

Shivani Goel  
Name

QC  
Title

Shivani Goel 08 Nov 2021  
Signature/Date