

Product Number:	101735
Lot Number:	502849
Expiration:	2-Mar-2022

Component	Item Number	Lot Number
SBG Dilution Buffer	100376	108322
SBG Concentrate	103397	106103

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 02Apr2021  
Signature/Date