

Product Number:	101735
Lot Number:	503193
Expiration:	19-Jul-2022

Component	Item Number	Lot Number
SBG Dilution Buffer	100376	123912
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 22Oct2021  
Signature/Date