

Product Number:	101735
Lot Number:	503087
Expiration:	13-Jul-2022

Component	Item Number	Lot Number
SBG Dilution Buffer	100376	120029
SBG Concentrate	103397	119416

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 04 Aug 2021
Signature/Date