

Product Number:	101735
Lot Number:	502486
Expiration:	3-Jun-2021

Component	Item Number	Lot Number
SBG Dilution Buffer	100376	015508
SBG Concentrate	103397	017607

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 24 Aug 2020
Signature/Date