

Product Number:	101735
Lot Number:	502968
Expiration:	15-Mar-2022

Component	Item Number	Lot Number
SBG Dilution Buffer	100376	112808
SBG Concentrate	103397	107425

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 03 Jun 2024
Signature/Date