

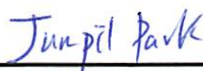


Product Number:	101735
Lot Number:	502987
Expiration:	24-Mar-2022

Component	Item Number	Lot Number
SBG Dilution Buffer	100376	112808
SBG Concentrate	103397	108324

Laboratory Analysis

Means of Analysis	Characteristic	Result
HD-1 Analyzer	Functional Test	PASS
SBG Concentrate Concentration (nM)		50

Review/Approval

Name Title Signature/Date