

Product Number:	101361
Lot Number:	504066
Expiration:	30-Oct-2024

Component	Item Number	Lot Number
SBG Dilution Buffer	100376	330321
SBG Concentrate	103397	330603
RGP Reagent	103159	N/A

¹ RGP is not Kit Lot Specific

Laboratory Analysis

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

Review/Approval

Sonia Khillan Senior QC Analyst Sonia Khillan 29 Nov 2023
Name Title Signature/Date