

Product Number:	101361
Lot Number:	503077
Expiration:	1-Jun-2022

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	120023	N/A
SBG Concentrate	103397	119416	50 nM
RGP Reagent	103159	N/A ¹	N/A

¹ RGP is not Kit Lot Specific

Review/Approval

Shivani Gael
Name

QC

Title

Shivani Gael 29 Jul 2021
Signature/Date