

Product Number:	101361
Lot Number:	502754
Expiration:	15-Dec-2021


Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	100820	N/A
SBG Concentrate	103397	101912	50 nM
RGP Reagent	103159	N/A ¹	N/A

¹ RGP is not Kit Lot Specific

Review/Approval

Junpil Park
Name

Supervisor, QC
Title

 23 Feb 2021
Signature/Date