

Product Number:	101361
Lot Number:	502617
Expiration:	27-Aug-2021

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	026623	N/A
SBG Concentrate	103397	024007	50 nM
RGP Reagent	103159	N/A ¹	N/A

¹ RGP is not Kit Lot Specific

Review/Approval

Shivani Goel
Name

Manager, QC
Title

Shivani Goel 17Nov2020
Signature/Date