

Product Number:	101361
Lot Number:	503432
Expiration:	26-Jan-2023

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	208905	N/A
SBG Concentrate	103397	204911	50 nM
RGP Reagent	103159	N/A ¹	N/A

¹ RGP is not Kit Lot Specific

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 04Apr2022

Name

Title

Signature/Date