

Product Number:	101361
Lot Number:	502912
Expiration:	15-Mar-2022

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	108322	N/A
SBG Concentrate	103397	107422	50 nM
RGP Reagent	103159	N/A ¹	N/A

¹ RGP is not Kit Lot Specific

Review/Approval

Shivani Goel
Name

QC

Title

Shivani Goel 04 May 2021
Signature/Date