

Product Number:	101361
Lot Number:	502789
Expiration:	11-Jan-2022

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	104727	N/A
SBG Concentrate	103397	101913	50 nM
RGP Reagent	103159	N/A ¹	N/A

¹ RGP is not Kit Lot Specific

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 10 Mar 2021
Signature/Date