

Product Number:	101361
Lot Number:	502352
Expiration:	27-Jan-2021

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	005223	N/A
SBG Concentrate	103397	002716	50 nM
RGP Reagent	103159	N/A ¹	N/A

¹ RGP is not Kit Lot Specific

Review/Approval

Shivam Goel
Name

QC Manager
Title

Shivam Goel 30 Apr 2020
Signature/Date