

Product Number:	101361
Lot Number:	502419
Expiration:	21-Feb-2021

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	005225	N/A
SBG Concentrate	103397	013908	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 09 Jun 2020
Signature/Date