

Product Number:	101361
Lot Number:	502528
Expiration:	03-Jun-2021

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	015511	N/A
SBG Concentrate	103397	017607	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 11 Sep 2020
Signature/Date