

Product Number:	101361
Lot Number:	502474
Expiration:	27-Jan-2021

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	015508	N/A
SBG Concentrate	103397	002715	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 20/1/2020
Signature/Date