

Product Number:	101361
Lot Number:	503012
Expiration:	24-Mar-2022

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	115221	N/A
SBG Concentrate	103397	108324	50 nM
RGP Reagent	103159	N/A ¹	N/A

¹ RGP is not Kit Lot Specific

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 24 Jun 2021
Signature/Date