

Product Number:	101361
Lot Number:	503439
Expiration:	19-Oct-2022

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	203113	N/A
SBG Concentrate	103397	129210	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 13Apr2022
Signature/Date