

Product Number:	101361
Lot Number:	502188
Expiration:	29-Oct-2020

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	930220	N/A
SBG Concentrate	103397	930316	50 nM
RGP Reagent	103159	N/A ¹	N/A

¹ RGP is not Kit Lot Specific

Review/Approval

Patrizia Stadler
Name

Assoc. Director, QC
Title

P. Stadler *09 Jan 2020*
Signature/Date