

Product Number:	101361
Lot Number:	502098
Expiration:	24-Jul-2020

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	921117	N/A
SBG Concentrate	103397	920511	50 nM
RGP Reagent	103159	N/A ¹	N/A

¹ RGP is not Kit Lot Specific

Review/Approval

Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i> 01 NOV 2019
Name	Title	Signature/Date