

Product Number:	101361
Lot Number:	503507
Expiration:	22-Feb-2023

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	211404	N/A
SBG Concentrate	103397	209717	50 nM
RGP Reagent	103159	N/A ¹	N/A

¹ RGP is not Kit Lot Specific

Review/Approval

Matt Sment
Name

QC
Title

MSA 02Jan2022
Signature/Date