

Product Number:	101361
Lot Number:	504190
Expiration:	25-Jul-2025

Component	Part Number	Lot Number	Control Number	Concentration
SBG Dilution Buffer	100376	402514	N/A	N/A
SBG Concentrate	103397	319211	319211A	50 nM
RGP Reagent	103159	N/A ¹	N/A	N/A

¹ Reagents are not Kit Lot Specific

Review/Approval

ROBERT SOMOGYI QC MANAGER

 22 MAR 2023

Name

Title

Signature/Date