

Product Number:	101361
Lot Number:	504505
Expiration:	15-Jul-2026

Component	Part Number	Lot Number	Control Lot Number	Concentration
SBG Dilution Buffer	100376	501523	501523A	N/A
SBG Concentrate	103397	502104	502104A	50 nM
RGP Reagent	103159	N/A ¹	N/A	N/A

¹ Reagents are not Kit Lot Specific

Review/Approval

William O
Name

Quality Systems Engineer I
Title

William O 06FEB2025
Signature/Date