

Product Number:	101361
Lot Number:	504460
Expiration:	3-Mar-2026

Component	Part Number	Lot Number	Control Lot Number	Concentration
SBG Dilution Buffer	100376	426913	N/A	N/A
SBG Concentrate	103397	319213	319213A	50 nM
RGP Reagent	103159	N/A ¹	N/A	N/A

¹ Reagents are not Kit Lot Specific

Review/Approval

TAM LEE DOC CONTROL II, QA Tam Lee 19 NOV 2024
 Name Title Signature/Date