

Product Number:	101361
Lot Number:	504340
Expiration:	28-Sep-2025

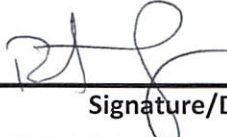
Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	417210	N/A
SBG Concentrate	103397	408804	50 nM
RGP Reagent	103159	N/A ¹	N/A

¹ Reagents are not Kit Lot Specific

Review/Approval

Robert Somogyi

QC Manager


05 AUG 2024

Name

Title

Signature/Date