

Product Number:	101361
Lot Number:	504174
Expiration:	17-Jul-2025

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	402514	N/A
SBG Concentrate	103397	401702	50 nM
RGP Reagent	103159	N/A ¹	N/A

¹ Reagents are not Kit Lot Specific

Review/Approval

ROBERT SOMOGYI	QC MANAGER	 07 MAR 2024
Name	Title	Signature/Date