

Product Number:	101361
Lot Number:	504360
Expiration:	20-Dec-2025

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	417210	N/A
SBG Concentrate	103397	420404	50 nM
RGP Reagent	103159	N/A <sup>1</sup>	N/A

<sup>1</sup> Reagents are not Kit Lot Specific

Review/Approval

*Dipesh Patel*  
Name

QA-IT  
Title

*[Signature]* / 23 Aug 2024  
Signature/Date