

<b>Product Number:</b>	101361
<b>Lot Number:</b>	504397
<b>Expiration:</b>	9-Jan-2026

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	424714	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

Dipak Patel      Doc Control - II QA

Name

Title

ant / 25 Sep 2024

Signature/Date