

<b>Product Number:</b>	101361
<b>Lot Number:</b>	502503
<b>Expiration:</b>	18-May-2021

Component	Part Number	Lot Number	Concentration
<b>SBG Dilution Buffer</b>	100376	015508	N/A
<b>SBG Concentrate</b>	103397	013909	50 nM
<b>RGP Reagent</b>	103159	N/A <sup>1</sup>	N/A

<sup>1</sup> RGP is not Kit Lot Specific

**Review/Approval**

Shivani Goel

QC Manager

*Shivani Goel 18 Aug 2020*

**Name**

**Title**

**Signature/Date**