

<b>Product Number:</b>	101735
<b>Lot Number:</b>	502402
<b>Expiration:</b>	2-May-2021

<b>Component</b>	<b>Item Number</b>	<b>Lot Number</b>
<b>SBG Dilution Buffer</b>	100376	012306
<b>SBG Concentrate</b>	103397	012511

<b>SBG Concentrate Concentration (nM)</b>	50
-----------------------------------------------	----

Review/Approval

Shivani Goel  
Name

QC Manager  
Title

*Shivani Goel* 02 Jun 2020  
Signature/Date