

<b>Product Number:</b>	101735
<b>Lot Number:</b>	501501
<b>Expiration:</b>	6-Aug-2019

<b>Component</b>	<b>Item Number</b>	<b>Lot Number</b>
<b>SBG Concentrate</b>	100439	821806
<b>SBG Dilution Buffer</b>	100376	829912

<b>SBG Concentrate Concentration (nM)</b>	144
---	-----

**Review/Approval**

Sivasankari Murugar      QC Supervisor  
\_\_\_\_\_  
Name                                      Title

*[Signature]*      19 NOV 2018  
\_\_\_\_\_  
Signature/Date

Statement: The above product was tested using Quanterix procedures and conforms to Quanterix's specifications.