

<b>Product Number:</b>	101735
<b>Lot Number:</b>	501399
<b>Expiration:</b>	10-Jul-2019

Component	Item Number	Lot Number
<b>SBG Concentrate</b>	100439	819109
<b>SBG Dilution Buffer</b>	100376	821410

<b>SBG Concentrate Concentration (nM)</b>	151
---	-----

**Review/Approval**

Sivasankari Murugar QC Supervisor  
Name Title

 05 SEP 2018  
Signature/Date

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