

<b>Product Number:</b>	101735
<b>Lot Number:</b>	501385
<b>Expiration:</b>	23-May-2019

Component	Item Number	Lot Number
SBG Concentrate	100439	817809
SBG Dilution Buffer	100376	MF-052318-12

SBG Concentrate Concentration (nM)	154
------------------------------------	-----

Review/Approval

Sivasankari Murugar QC Supervisor  
Name Title

  
Signature/Date

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