

<b>Product Number:</b>	101735
<b>Lot Number:</b>	501226
<b>Expiration:</b>	2-Apr-2019

<b>Component</b>	<b>Item Number</b>	<b>Lot Number</b>
SBG Dilution Buffer	100376	MF-041518-05
SBG Concentrate	100439	80925

SBG Concentrate Concentration (nM)	144
---------------------------------------	-----

Review/Approval

Sivasankari Murugar QC Supervisor  
Name Title

  
Signature/Date 26 APR 2018

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