

<b>Product Number:</b>	101735
<b>Lot Number:</b>	501587
<b>Expiration:</b>	29-Aug-2019

Component	Item Number	Lot Number
SBG Concentrate	100439	824116
SBG Dilution Buffer	100376	834903

SBG Concentrate Concentration (nM)	155
---------------------------------------	-----

Review/Approval	Patrizia Stadler Name	Associate Director, QC Title	P. Stadler Signature/Date	24 Jan 2019 Date
-----------------	--------------------------	------------------------------------	------------------------------	---------------------

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