

Product Number:	101735
Lot Number:	501717
Expiration:	22-Jan-2020

Component	Item Number	Lot Number
SBG Concentrate	103397	907013
SBG Dilution Buffer	100376	902215

SBG Concentrate Concentration (nM)	50
---------------------------------------	----

Review/Approval	Patrizia Stadler	Associate Director, QC	<i>P. Stadler</i>	<i>18 Mar 2019</i>
	Name	Title	Signature/Date	

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