

Product Number:	101735
Lot Number:	501645
Expiration:	19-Nov-2019

Component	Item Number	Lot Number
SBG Concentrate	103397	832316
SBG Dilution Buffer	100376	834903

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

Review/Approval	Patrizia Stadler	Associate Director, QC	<i>P. Stadler</i>	07 Feb 2019
	Name	Title	Signature/Date	

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