

<b>Product Number:</b>	101735
<b>Lot Number:</b>	501688
<b>Expiration:</b>	22-Jan-2020

Component	Item Number	Lot Number
SBG Concentrate	103397	902932
SBG Dilution Buffer	100376	902215

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

Review/Approval	Patrizia Stadler Name	Associate Director, QC Title	<i>p. Stodler</i> <i>1 March 2019</i> Signature/Date
-----------------	--------------------------	------------------------------------	---

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