

<b>Product Number:</b>	101735
<b>Lot Number:</b>	501746
<b>Expiration:</b>	11-Mar-2020

Component	Item Number	Lot Number
SBG Concentrate	103397	907016
SBG Dilution Buffer	100376	907014

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

Review/Approval	Patrizia Stadler Name	Associate Director, QC Title	<i>P. Stadler</i> 29 Mar 2019 Signature/Date
-----------------	--------------------------	------------------------------------	---

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