

Product Number:	103397
Lot Number:	925305
Expiration:	10-Sep-2020
Storage:	2 - 8°C

SBG Concentrate Concentration (nM)	141
---------------------------------------	-----

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

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