

Product Number:	101735
Lot Number:	502054
Expiration:	30-Jul-2020

Component	Item Number	Lot Number
SBG Dilution Buffer	100376	921119
SBG Concentrate	103397	925509

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

Review/Approval

Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i>	04 OCT 2019
Name	Title	Signature/Date	