

Product Number:	101735
Lot Number:	502175
Expiration:	29-Oct-2020

Component	Item Number	Lot Number
SBG Dilution Buffer	100376	930220
SBG Concentrate	103397	930316

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

Review/Approval

Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i>	06 Jan 2020
Name	Title	Signature/Date	