

Product Number:	101735
Lot Number:	501877
Expiration:	15-Apr-2020

Component	Item Number	Lot Number
SBG Dilution Buffer	100376	916412
SBG Concentrate	103397	910511

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

Review/Approval

Patrizia Stadler	Assoc. Director, QC	<i>p. stadler</i>	27 JUN 2019
Name	Title	Signature/Date	