

<b>Product Number:</b>	101735
<b>Lot Number:</b>	502118
<b>Expiration:</b>	12-Sep-2020

Component	Item Number	Lot Number
<b>SBG Dilution Buffer</b>	100376	930219
<b>SBG Concentrate</b>	103397	925509

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

Review/Approval

Patrizia Stadler

Assoc. Director, QC

*P. Stadler*

14 Nov 2019

Name

Title

Signature/Date