

<b>Product Number:</b>	101735
<b>Lot Number:</b>	502096
<b>Expiration:</b>	12-Sep-2020

Component	Item Number	Lot Number
SBG Dilution Buffer	100376	926017
SBG Concentrate	103397	925509

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

Review/Approval

Patrizia Stadler  
Name

Assoc. Director, QC  
Title

*p.stadler*

29 Oct 2019  
Signature/Date