

<b>Product Number:</b>	101735
<b>Lot Number:</b>	502211
<b>Expiration:</b>	21-Nov-2020

<b>Component</b>	<b>Item Number</b>	<b>Lot Number</b>
<b>SBG Dilution Buffer</b>	100376	932513
<b>SBG Concentrate</b>	103397	935218

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

Review/Approval      Patrizia Stadler      Assoc. Director, QC      P. Stadler      29 Jan 2020  
Name      Title      Signature/Date