

Product Number:	101361
Lot Number:	503460
Expiration:	8-Feb-2023

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	208911	N/A
SBG Concentrate	103397	204912	50 nM
RGP Reagent	103159	N/A ¹	N/A

¹ RGP is not Kit Lot Specific

Review/Approval

<p><i>Richa Shrestha</i> <u>Shivani Goel</u></p>	<p><i>QC Supervisor</i> <u>QC Manager</u></p>	<p><i>[Signature]</i> <u>Signature/Date</u></p>
<p>Name <i>w/22Apr22</i></p>	<p>Title <i>w/22Apr22</i></p>	<p><i>22Apr 2022</i></p>