

|                        |             |
|------------------------|-------------|
| <b>Product Number:</b> | 101361      |
| <b>Lot Number:</b>     | 502871      |
| <b>Expiration:</b>     | 11-Feb-2022 |

| Component           | Part Number | Lot Number       | Concentration |
|---------------------|-------------|------------------|---------------|
| SBG Dilution Buffer | 100376      | 108322           | N/A           |
| SBG Concentrate     | 103397      | 104730           | 50 nM         |
| RGP Reagent         | 103159      | N/A <sup>1</sup> | N/A           |

<sup>1</sup> RGP is not Kit Lot Specific

Review/Approval

*Junpil Park*  
Name

QC Supervisor  
Title

*[Signature]* 14 Apr 2021  
Signature/Date