

<b>Product Number:</b>	101735
<b>Lot Number:</b>	502532
<b>Expiration:</b>	3-Jun-2021

Component	Item Number	Lot Number
SBG Dilution Buffer	100376	015511
SBG Concentrate	103397	017607

**Laboratory Analysis**

Means of Analysis	Characteristic	Result
HD-1 Analyzer	Functional Test	PASS
SBG Concentrate Concentration (nM)		50

Review/Approval

Shivani Goel  
Name

QC Manager  
Title

Shivani Goel 11 Sep 2020  
Signature/Date