

Product Number:	101735
Lot Number:	503141
Expiration:	28-Jun-2022

Component	Item Number	Lot Number
SBG Dilution Buffer	100376	121719
SBG Concentrate	103397	120031

**Laboratory Analysis**

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

Review/Approval

*Shivani Goel*

Name

QC

Title

*Shivani Goel 15 Sep 2021*

Signature/Date