

Product Number:	103650
Lot Number:	117921
Expiration:	28-Sep-2021
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
HD-1 Analyzer	PASS

SBG Concentrate Concentration (nM)	152
---------------------------------------	-----

Review/Approval

Shivani Goel

Name

QC

Title

Shivani Goel 19 Jul 2021

Signature/Date