

Product Number:	103650
Lot Number:	034212
Expiration:	7-Mar-2021
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
HD-1 Analyzer	PASS

SBG Concentrate Concentration (nM)	143
---------------------------------------	-----

Review/Approval

Junpil Park
Name

Supervisor, QC
Title

 08 Jan 2021
Signature/Date