

Product Number:	103397
Lot Number:	225614
Expiration:	13-Sep-2023
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
HD-1 Analyzer	PASS

SBG Concentrate Concentration (nM)	50
------------------------------------	----

Review/Approval Sonia Khillan QC Manager Sonia Khillan / 18 JAN 2023
Name Title Signature/Date