

Product Number:	103397
Lot Number:	308712
Expiration:	28-Mar-2024
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
HD-1 Analyzer	PASS

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

Review/Approval

Sonia K. Miller
Name

QC Manager
Title

Sonia K. Miller 20 JUN 2023
Signature/Date