

Product Number:	103397
Lot Number:	408804
Expiration:	28-Sep-2025
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	PASS / FAIL
Functionality	HD(X/1) Analyzer	PASS

RESULT

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

Review/Approval

ROBERT SOMOGYI

QC MANAGER

[Signature] 05 APR 2024

Name

Title

Signature/Date