

Product Number:	103397
Lot Number:	315712
Control Number:	315712A
Expiration:	6-Jun-2026
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Spectrophotometer	PASS

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

Review/Approval

ROBERT SAMOGRY

Name

QC MANAGER

Title

[Signature] 07 AUG 2023

Signature/Date