

Product Number:	104360
Lot Number:	315806
Control Lot Number:	315806A
Expiration:	7-Jun-2029
Storage:	-80°C

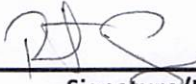
Laboratory Analysis

Characteristic	Result
Concentration	1 mg/ml

Review/Approval

Robert Somogyi

QC Manager

 05 MAR 2025

Name

Title

Signature/Date