

Product Number:	102243
Lot Number:	321204
Expiration:	31-Jul-2025
Storage:	-80°C

Laboratory Analysis

Means of Analysis	Characteristic	(mg/mL)
SR-X Analyzer	Nominal Concentration	910.4

Review/Approval

ROBERT SOMOGYI
Name

QC MANAGER
Title

 18 JAN 2024
Signature/Date