

Product Number:	102619
Lot Number:	329214
Expiration:	19-Oct-2025
Storage:	- 20°C

Laboratory Analysis

Concentration (pg/ml) (Tested with N2PB)	12275
Concentration (pg/ml) (Tested with N4PE)	15619

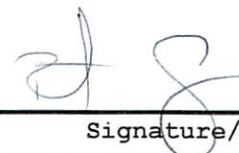
Review/Approval

ROBERT SOMOGYI

Name

QC MANAGER

Title

 23 JAN 2024

Signature/Date