

Product Name:	Neuro 4-Plex E Control 0
Product Number:	103846
Lot Number:	114605
Paired Reagent Kit Lot Number:	502959
Expiration:	26-May-2023
Platform(s):	HD-X
Storage Condition:	-80°C

Laboratory Analysis			
Characteristic	Mean concentration Range		
Means of Analysis	HD-X Analyzer		
Result	Plex	Mean Range (pg/mL)	Ranges shown are generated for internal use only. Customer should generate their own control ranges.
	Abeta 40	5.24-9.55	
	Abeta 42	2.07-3.10	
	GFAP®	102-153	
	NF-light®	12.2-18.7	

Review/Approval

Junpil Park

QC Supervisor

 04 Jun 2024

Name

Title

Signature/Date