

Product Name:	Neuro 4-Plex E Control 0
Product Number:	103846
Lot Number:	028115
Paired Reagent Kit Lot Number:	502542
Expiration:	7-Oct-2022
Platform(s):	HD-X & SR-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	6.12-9.17	
	Abeta 42	2.19-3.28	
	GFAP®	86.0-129	
	NF-light®	9.05-17.6	

Review/Approval

Patrizia Stadler Associate Director, QC *P. Stadler* 22Oct2020
 Name Title Signature/Date