

Neuro 4-Plex E Control 0	
Product Number:	103846
Lot Number:	206909
Paired Reagent Kit : Neuro 4-Plex E Bulk Kit (item 103971)	
Lot Number:	503415
Expiration:	24-Nov-2022
Platform(s):	HD-X
Storage Condition:	-80°C



*Data below represents
results generated on
the Simoa® HD-X
Analyzer.*

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	5.26-10.0	
	Abeta 42	1.75-2.62	
	GFAP®	64.6-96.9	
	NF-light®	7.10-10.7	

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 25 Mar 2022

Name

Title

Signature/Date