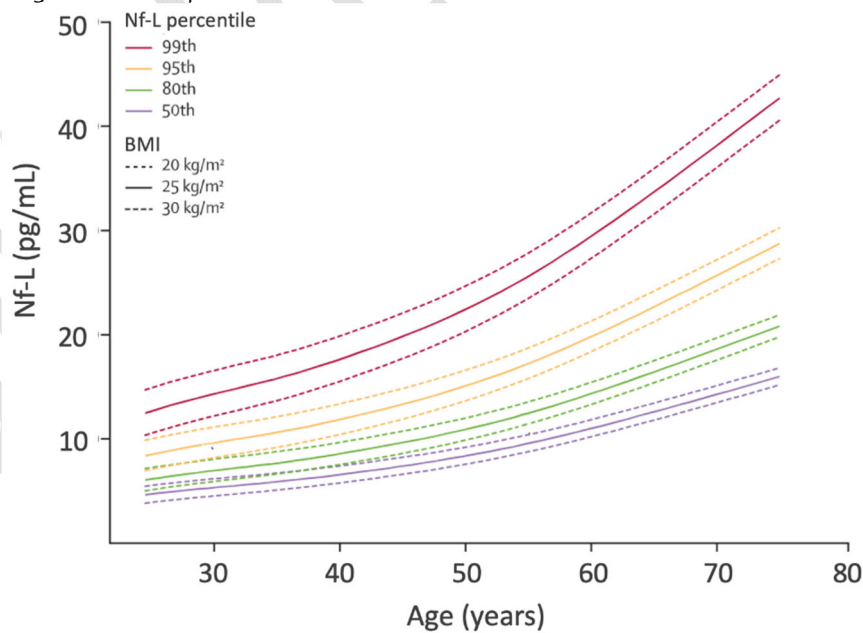


Patient Name		
Sex	Age	Date of Birth <small>DD-MMM-YYYY</small>
Sample Number		Sample Type <input type="checkbox"/> Serum
Provider Name		Collection Date <small>DD-MMM-YYYY</small>
Organization		Collection Time <small>HH:MM</small>
Secure Fax#:		Date Received <small>DD-MMM-YYYY</small>
		Report Date <small>DD-MMM-YYYY</small>

	Result (pg/mL)		
104291 Quanterix Nf-L Test, Serum	XX.XX		
Interpretation: Nf-L elevation in blood can be due to many different causes, including neurodegenerative diseases or a head impact. Nf-L results should only be used in conjunction with other clinical information when evaluating patients. Various demographic, lifestyle, and comorbidity factors can influence Nf-L levels in serum. The largest demographic variable is age, necessitating age-adjusted reference intervals. A reference range study was conducted across 113 presumed normal healthy individuals varying in age from 19 to 69 years old. Table 1 exhibits median values and expected ranges of serum Nf-L in different age groups based on 5th percentile through 95th percentiles as estimated from Anderson-Darling fitting.	Age	Median, pg/ml	Range, pg/mL
	19-39	4.86	2.47-7.25
	40-49	7.01	4.44-9.58
	50-59	11.02	4.79-17.25
	60-69	13.12	3.97-22.27

Nf-L test results can be further visualized on the chart below which is based on a normative study of over 10,000 samples and exhibits both the age and BMI dependence of serum Nf-L.¹



Test Information:

The Quanterix Nf-L Test is a digital immunoassay for the quantitative determination of neurofilament light chain (Nf-L) in human serum. Nf-L is a 68 kDa cytoskeletal intermediate filament protein that is expressed in neurons. It associates with the 125 kDa neurofilament medium and the 200 kDa neurofilament heavy chains to form neurofilaments.² Neurofilaments are major components of the neuronal cytoskeleton and are believed to function primarily to provide structural support for the axon and to regulate axon diameter.³ Neurofilaments can be released in significant quantities into the cerebrospinal fluid (CSF) following axonal damage or neuronal degeneration. A fraction of these proteins diffuses into the blood, where concentrations are typically 50 to 100-fold lower than in CSF.⁴ Nf-L elevations have been demonstrated with traumatic brain injury, multiple sclerosis, frontotemporal dementia, Alzheimer's disease, and other neurodegenerative diseases.⁵⁻⁸

REFERENCES

1. Benkert P, Meier S, Schaedelin S, et al. Serum neurofilament light chain for individual prognostication of disease activity in people with multiple sclerosis: a retrospective modelling and validation study. *Lancet Neurol.* 2022;21(3):246-257.
2. Yuan A, Rao MV, Veeranna, Nixon RA. Neurofilaments at a glance. *J Cell Sci.* 2012;125(Pt 14):3257-3263. doi:10.1242/jcs.104729
3. Lalonde R, Strazielle C, et al. Neurobehavioral characteristics of mice with modified intermediate filament genes. *Rev Neurosci* 2003; 14 (4): 369–85.
4. Gaiottino J, Norgren N, Dobson R, et al. Increased neurofilament light chain blood levels in neurodegenerative neurological diseases. *PLoS One.* 2013;8(9):e75091. Published 2013 Sep 20.
5. Shahim P, Gren M, Liman V, Andreasson U, Norgren N, Tegner Y, Mattsson N, Andreassen N, Öst M, Zetterberg H, Nellgård B, Blennow K. Serum neurofilament light protein predicts clinical outcome in traumatic brain injury. *Sci Rep.* 2016 Nov 7;6:36791.
6. Disanto G, Barro C, Benkert P, et al. Serum Neurofilament light: A biomarker of neuronal damage in multiple sclerosis. *Ann Neurol.* 2017;81(6):857-870.
7. Meeter LH, Dopfer EG, et al. Neurofilament light chain: a biomarker for genetic frontotemporal dementia. *Ann Clin Transl Neurol* 2016;3(8):623#36.
8. Giacomucci G, Mazzeo S, Bagnoli S, Ingannato A, Leccese D, Berti V, Padiglioni S, Galdo G, Ferrari C, Sorbi S, Bessi V, Nacmias B. Plasma neurofilament light chain as a biomarker of Alzheimer's disease in subjective cognitive decline and mild cognitive impairment. *J Neurol.* 2022 Aug;269(8):4270-4280.

Quanterix Laboratory Director: *Timothy Skelton, M.D., Ph.D., ABPD*