

NF-light® (Neurofilament light)

Quanterix™

UmanDiagnostics
A Quanterix Company

NF-light® ELISA CE-IVD

Intended use:

NF-light® ELISA is an invitro diagnostic device intended for quantitative determinations of human Neurofilament light (NF-L) protein in cerebrospinal fluid (CSF). The result is used to aid the diagnosis of neurological diseases such as amyotrophic lateral sclerosis (ALS), multiple sclerosis (MS), dementias and Parkinson's (PD). The kit is intended for professional use.

In addition, the NF-light® ELISA can be used for research using samples containing NF-L from rat, mouse, bovine and macaques sources as the antibodies in the assay recognizes NF-L from these species as well.

Technology: 2-site solid phase sandwich ELISA

Contents: Materials sufficient for 96 determinations including the Standard curve

Assay procedure:

- Dilute the CSF samples (50µl) with equal amount (1+1) of Sample diluent
- Wash the wells to be used with Wash buffer (3x300 µL).
- Add 100 µL of each Standard and sample in duplicate. Incubate 1 hour at RT with agitation
- Wash the wells with Wash buffer
- Add 100 µL of freshly diluted Tracer antibody to each well. Incubate 45 minutes at RT with agitation
- Wash the wells with Wash buffer
- Add 100 µL of newly diluted Conjugate to each well. Incubate 30 minutes at RT with agitation
- Wash the wells with Wash buffer
- Add 100 µL of TMB to each well. Incubate 15 minutes at RT with agitation
- Add 50 µL of Stop reagent to each well and read the absorbance at 450 nm (reference wavelength 620-650 nm)

Assay performance of the CE-certified NF-light® ELISA

| | |
|-------------------------|---|
| Detection limit: | 33ng/L |
| Measuring range: | 100 to 10 000 ng/L |
| Precision: | Intra-assay CV < 5% Inter-assay CV < 10% |
| Hook effect: | No Hook effect below 10 000 ng/L |

| | | |
|--------------------|----|--|
| References: | 1) | Rosengren et al. J Neurochem. 1996. 67(5):2013-8 |
| | 2) | Norgren et al. Brain Res. 2003. 987(1):25-31 |
| | 3) | Norgren et al. Neurology. 2004. 63(9):1586-90 |
| | 4) | Norgren et al. Brain Res Bull. 2005. 67(4):264-8 |



Contact information:

UmanDiagnostics AB, tel. no: +46 90 777 880, Tvistevagen 48C, 90736 Umea, Sweden
e-mail: info@umandiagnosics.se, www.umandiagnosics.com