

Neopterin ELISA.

IMMUNOLOGY / CYTOKINES

Manufactured by

IBL International GmbH
Flughafenstrasse 52a
22335 Hamburg
Germany

Phone: +49 (0)40-53 28 91-0
Fax: +49 (0)40-53 28 91-11
Email: IBL@Tecan.com
www.tecan.com/immunoassays

Australia +61 3 9647 4100 **Austria** +43 62 46 89 330 **Belgium** +32 15 42 13 19 **China** +86 21 220 63 206 **France** +33 4 72 76 04 80 **Germany** +49 79 51 94 170
Italy +39 02 92 44 790 **Japan** +81 44 556 73 11 **Netherlands** +31 18 34 48 17 4 **Nordic** +46 8 750 39 40 **Singapore** +65 644 41 886 **Spain** +34 93 93 595 25 31
Switzerland +41 44 922 89 22 **UK** +44 118 9300 300 **USA** +1 919 361 5200 **Other countries** +41 44 922 81 11

Tecan Group Ltd. makes every effort to include accurate and up-to-date information; however, it is possible that omissions or errors might have occurred. Tecan Group Ltd. cannot, therefore, make any representations or warranties, expressed or implied, as to the accuracy or completeness of the information provided. Changes can be made at any time without notice. All mentioned trademarks are protected by law. In general, the trademarks and designs referenced herein are trademarks, or registered trademarks, of Tecan Group Ltd., Maennedorf, Switzerland. A complete list may be found at www.tecan.com/trademarks. Product names and company names that are not contained in the list but are noted herein may be the trademarks of their respective owners. For technical details and detailed procedures of the specifications provided please contact your Tecan representative. This may contain reference to applications and products which are not available in all markets. Please check with your local sales representative.

© 2020 Tecan Trading AG, Switzerland, all rights reserved. For disclaimer and trademarks please visit www.tecan.com

www.tecan.com

 **TECAN.**

Neopterin EN 2021 V01

A general marker for cellular immune system activation.

An increase in Neopterin concentration marks the onset of an infection in the body before antibodies are present. As soon as the immune system is triggered by interferon- γ , Neopterin is released by macrophages (Figure 1).

PRODUCT

RE59321
Neopterin ELISA (CE, IVD)

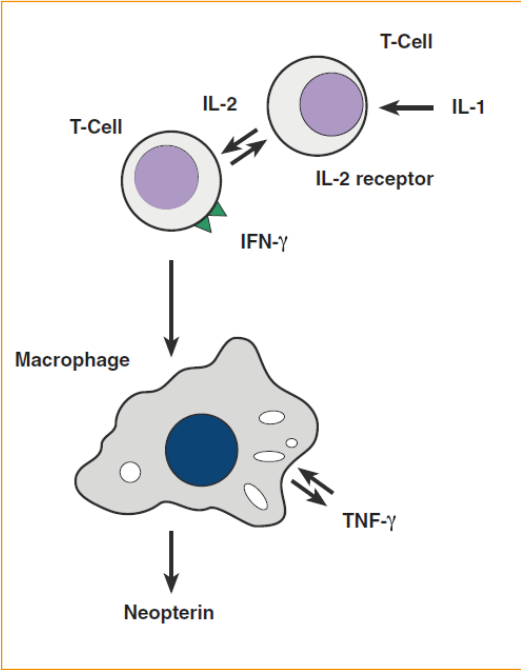


Figure 1: Release of Neopterin during cellular immune reaction.¹

EARLY DETECTION OF CELL IMMUNE REACTION

- Elevated Neopterin can be measured and is useful as¹:
- Aid for early diagnosis of viral infections (e.g. **SARS-CoV-2**, HIV.)
 - Differential diagnosis between viral and bacterial infections
 - Autoimmune diseases (Sjögren-syndrome, rheumatic diseases)
 - Early indication of complications in allograft rejection episodes
 - Follow-up control of chronic infections, therapy and treatment

The IBL International Neopterin ELISA is calibrated against the HPLC reference method. This calibration against external standards gives our products an extraordinarily high accuracy and precision from lot to lot.

Insights from Neopterin about the pathology of SARS-CoV-2
Synthesized upon stimulation of macrophages and dendritic cells, Neopterin is an early, well-characterized biochemical marker of an ongoing immune reaction to viral infection (Figure 2).²

There is evidence that Neopterin levels correlate with severity of COVID-19 progression and may be prognostic in COVID-19 patients.³⁻⁴ The Neopterin ELISA kit sensitively detects Neopterin prior to clinical symptoms of infection and thus, may be a tool to support an early recognition of SARS-CoV-2 infection.

PROGNOSIS, PROGRESSION AND THERAPY CONTROL

- Applications of Neopterin testing:
- Potential prognostic value in case of a COVID-19 infection
 - Aid for the detection of viral infections before onset of antibodies e.g. IgG, IgM
 - Monitoring the inflammation process
 - Monitoring the efficacy of an immune stimulatory therapy

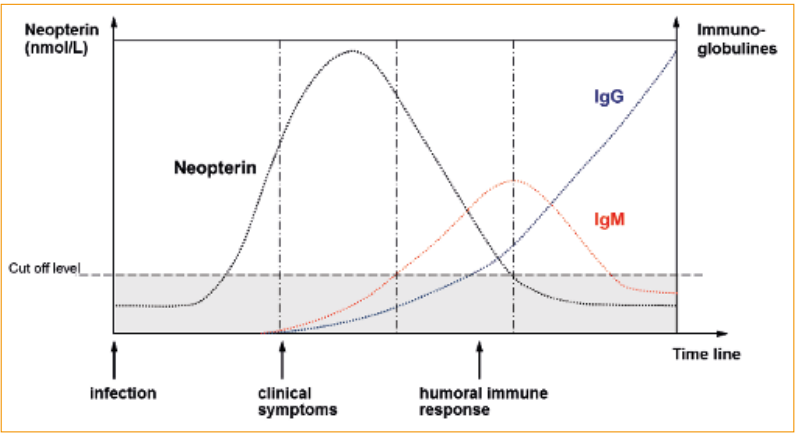


Figure 2: Schematic time course of Neopterin concentration in blood in relation to the rise and fall of antibody presence in response to viral infection.²

EASY TO PERFORM AND ACCURATE ELISA

The Neopterin ELISA is easy to perform, manually or on open ELISA processors, e.g. DSX, Etimax, EVOLyzer® and ThunderBolt® (The combined use of assays, process script and instrument has to be validated individually on site by each laboratory.). Expected values of Neopterin in serum are depicted in Table 1.

Serum	
nmol/L	ng/mL
> 10	< 2.5

Table 1: Normal values in serum measured in an apparently healthy population.

VERY GOOD CORRELATION TO HPLC

- Key benefits:
- Very good correlation to HPLC ($R^2=0.9974$)
 - Results within 2h
 - Validated for serum, plasma (EDTA) and urine. Further bodily fluids for research applications

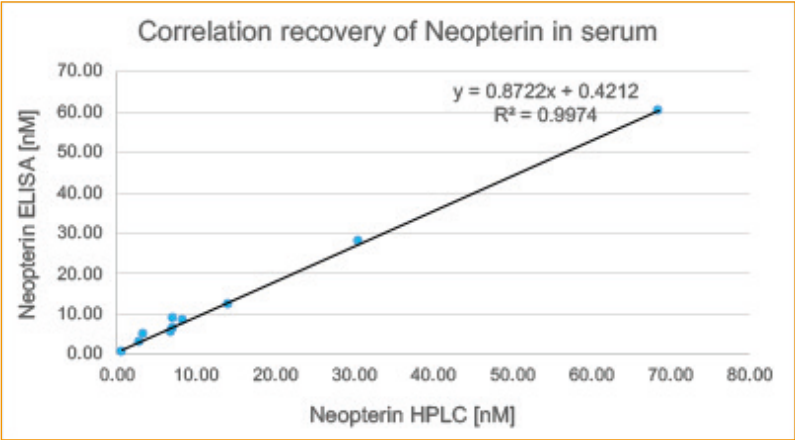


Figure 3: Method comparison between HPLC and IBL Neopterin ELISA (n=10, serum samples).

References:

1. Murr, C., Widner, B., Wirleitner, B., & Fuchs, D. (2002). Neopterin as a marker for immune system activation. Current drug metabolism, 3(2), 175-187.
2. Schennach, H., et al. (1994) Increased prevalence of IgM antibodies to Epstein-Barr virus and parvovirus B19 in blood donations with above-normal neopterin. Clinical chemistry 40: 2104. doi: 10.1128/CVI.00380-08
3. Robertson, J., et al. (2020) Serum neopterin levels in relation to mild and severe COVID-19. medRxiv. doi: 10.1101/2020.08.19.20178178
4. Ozger, H. S., et al. (2020). The prognostic role of neopterin in COVID-19 patients. Journal of Medical Virology. Doi: 10.1002/jmv.26472
5. Edén, A., et al. (2020) CSF biomarkers in patients with COVID-19 and neurological symptoms: A case series. Neurology. doi:10.1212/WNL.0000000000010977

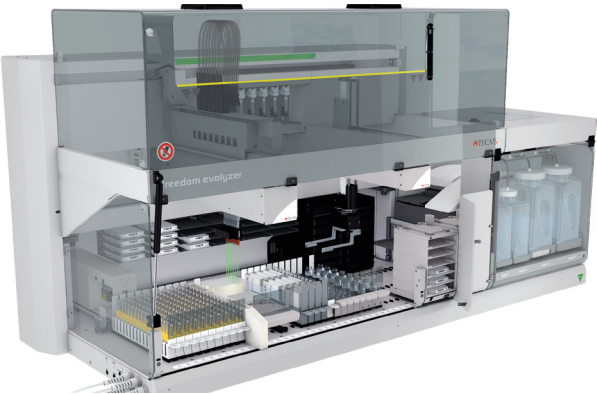


Figure 4: EVOLyzer®-Effective ELISA automation from Tecan