

# Nephrolyx



KDIGO 2024 CKD Guideline Gold Standard

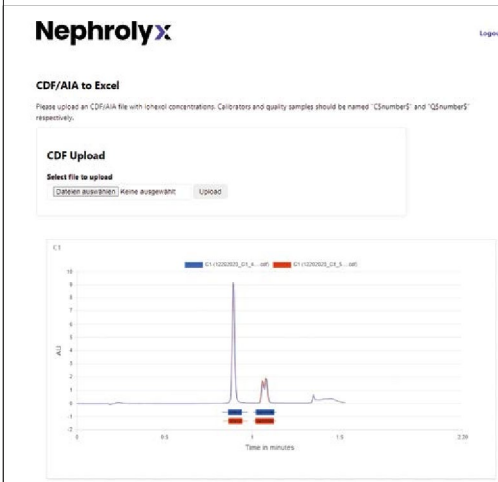
**True GFR Measurement  
by Iohexol Plasma Clearance**

# The Nephrolyx DNPx software platform enables smooth workflow integration of GFR measurement

## Digital Nephrology Platform (DNPx)

The GFR is determined by tracking the clearance of Iohexol by the patient's kidneys over time. Utilizing data from Iohexol administration, sampling times, and available patient-specific parameters, the mGFR is accurately calculated using a sophisticated algorithmic software application. The Nephrolyx DNPx offers flexible deployment options, including on-premises hosting or as a cloud solution within the facility's private network.

### Chromatogram



**Nephrolyx** Logout

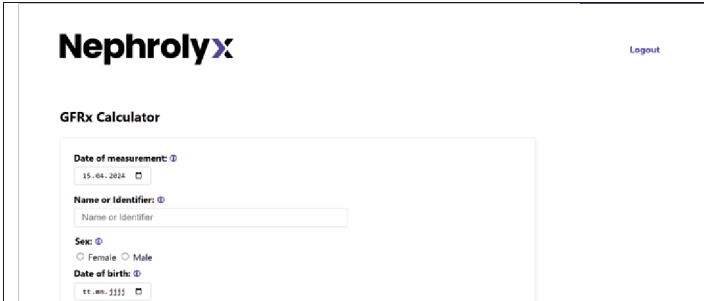
**CDF/AIA to Excel**  
Please upload an CDF/AIA file with Iohexol concentrations. Calibrators and quality samples should be named "C[number]" and "Q[number]" respectively.

**CDF Upload**  
Select file to upload

**Chromatogram**  
The plot shows Iohexol concentration (AU) on the y-axis (ranging from -2 to 10) versus Time in minutes on the x-axis (ranging from 0 to 3.00). Two peaks are visible, one at approximately 0.8 minutes and another at approximately 1.2 minutes.

UHPLC Settings	
<b>Solvent A</b>	UHPLC-grade water with 0.1 % formic acid
<b>Solvent B</b>	UHPLC-grade acetonitrile with 0.1 % formic acid
<b>Autosampler wash solution</b>	UHPLC-grade water with 10.0 % acetonitrile
<b>Flow</b>	1.0 ml/min
<b>Column temp.</b>	50 °C
<b>Inject volume</b>	10 µl
<b>Autosampler temperature</b>	4 °C
<b>Wavelength</b>	250 nm

### Patient/sample information input



**Nephrolyx** Logout

**GFRx Calculator**

**Date of measurement:** 15.04.2024

**Name or Identifier:**

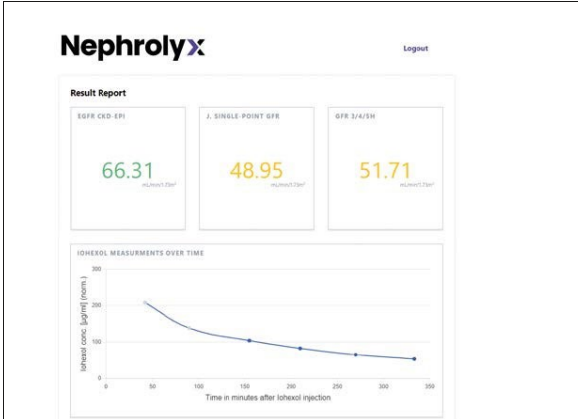
**Sex:**  Female  Male

**Date of birth:** 11.01.1951

Data transfer from LIMS and UHPLC instrument software can be automated & integrated.

- Manual input possible

### Result display



**Nephrolyx** Logout

**Result Report**

GFR CKD-EPI

66.31

mL/min/1.73m<sup>2</sup>

J. SINGLE POINT GFR

48.95

mL/min/1.73m<sup>2</sup>

GFR 3/4/5H

51.71

mL/min/1.73m<sup>2</sup>

**IOHEXOL MEASUREMENTS OVER TIME**

The graph shows Iohexol concentration (µg/ml (norm.)) on the y-axis (0 to 300) versus Time in minutes after Iohexol injection on the x-axis (0 to 350). The concentration starts at approximately 200 µg/ml at 0 minutes and decreases over time, reaching about 50 µg/ml at 350 minutes.

mGFR automatically calculated and reported:

- Area under the curve (AUC)
- Single and multiple-sample protocol GFR results
- Absolut or body surface indexed

# The Nephrolyx RUOx is a ready-to-use solution for direct kidney function measurement – mGFR

The test is intended for the in-vitro diagnostic use to quantify the serum iohexol concentration followed by the calculation of the resulting glomerular filtration rate (GFR) as an aid in the diagnosis, monitoring, and treatment of kidney diseases. The principle of the test is based on quantification using UHPLC with UV light detection. The chromatographic separation and elution of iohexol is performed using a linear gradient on a reversed-phase chromatography column.



## *Nephrolyx RUOx Test kit:*

*The kit is ready to use and allows up to 152 patient sample measurements in addition to calibration and QC measurements. It has an open-vial (in-use) stability of 14 days. Time to result: less than 3 hours; 90 seconds gradient.*

Nephrolyx RUOx test kit contains 6 calibrants and 3 quality control levels for use in the UHPLC analysis. It uses European Pharmacopeia Iohexol Certified Reference Standard as a primary reference material to establish a calibration hierarchy for both the end-user RUO calibrators and the quality control materials according to EN ISO 17511:2021.

## Key features and advantages

### **Providing confidence to physicians & patients**

Measuring Glomerular Filtration Rate (mGFR) through plasma clearance of the exogenous marker Iohexol is recognized as the gold standard by the latest Kidney Disease Improving Global Outcomes (KDIGO) 2024 CKD guideline. Consequently, Nephrolyx RUOx delivers the highest accuracy and precision in GFR measurement, enabling physicians to utilize precision medicine to enhance patient clinical outcomes.

### **Accuracy & ease of use**

The Nephrolyx RUOx provides direct GFR measurement with superior accuracy and precision over traditional biomarker-based estimates like serum creatinine, Cystatin C, and others. As a RUO product, it not only meets the highest regulatory standards but is also more user-friendly and easier to perform than lab-developed tests (LDTs) or nuclear medicine solutions. The kit is ready to use and allows up to 152 sample measurements in addition to calibration and QC measurements. It features an open-vial (in-use) stability of 14 days.

### **Rapid**

The Nephrolyx RUOx offers a streamlined workflow that delivers reliable results in less than three hours, making it faster than other mGFR methods.

## Reliable data

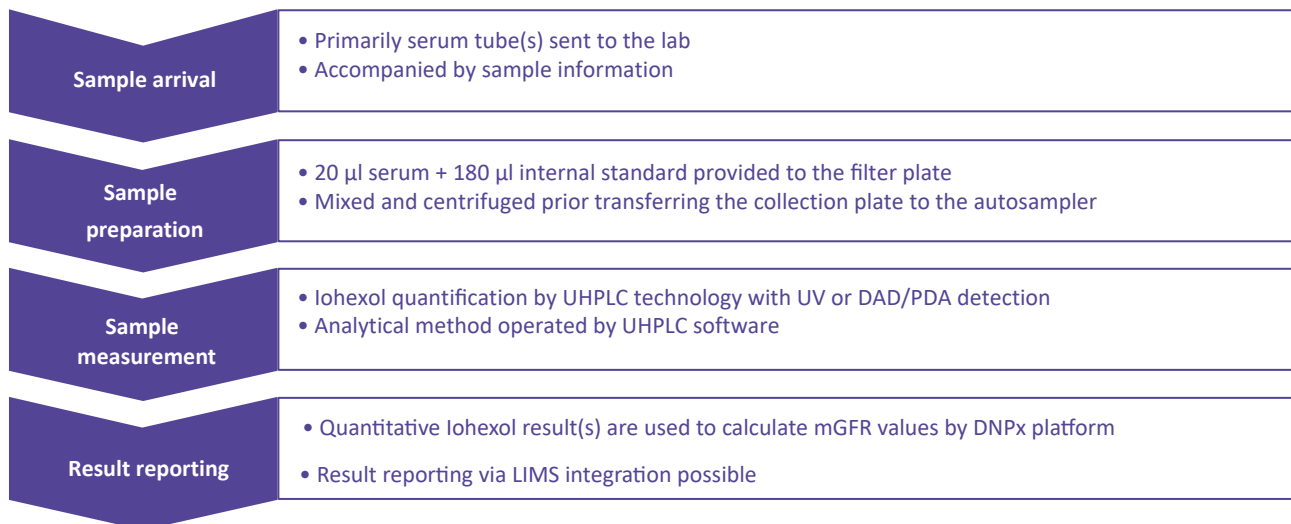
Operating Nephrolyx RUOx with a compatible UHPLC system enables excellent analytical performance in a fully regulatory compliant workflow (EU only).

## Lab perspective: From requirements to analytical workflow

Required UHPLC specifications	Required equipment	
<ul style="list-style-type: none"><li>• UPLC class system &gt; 500 bar binary pump system (e.g., Shimadzu, Hitachi, Waters, Agilent, Thermo Scientific)</li><li>• Autosampler with 4C colling function</li><li>• UV or DAD/PDA detector</li><li>• Column heater 50C</li><li>• 40-100 uL sample loop</li></ul>	<b>Analytical column</b>	PN 001310 "Nephrolyx UHPLC Column"
	<b>Guard column</b>	PN 001320 "Nephrolyx UHPLC Guard Column"
	<b>Platform shaker</b>	>= 1,000 rpm
	<b>Plate centrifuge</b>	= 1,500 G
	<b>Pipette for 20 µl</b>	Systematic error <= 1.0 %, random error <= 0.3 % With reverse pipetting support
	<b>Pipette for 180 µl</b>	Systematic error <= 0.8 %, random error <= 0.45 % With reverse pipetting support

The proprietary Nephrolyx guard and analytical columns are used to perform the test with UHPLC; they are not parts of the kit and need to be ordered additionally.

## Laboratory workflow



## Analytical specifications

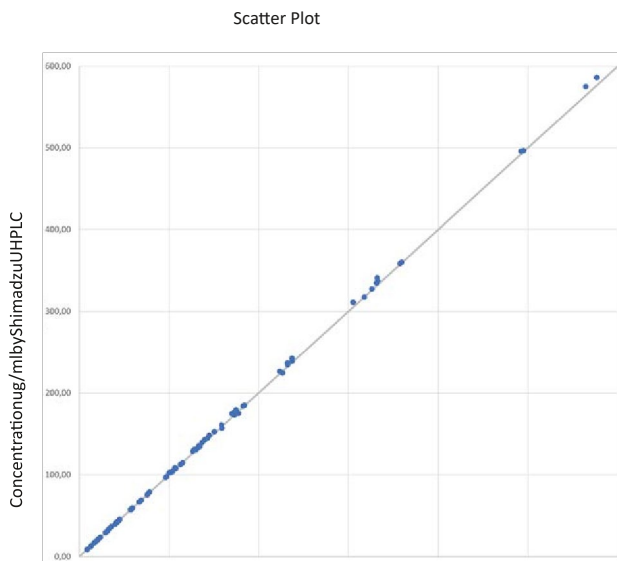


Figure 1 : Correlation between reference and Shimadzu Nexera X3

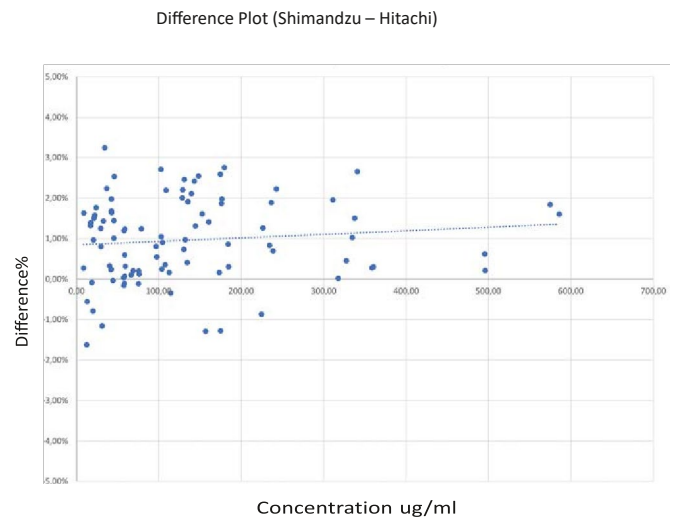


Figure 2 : Difference plot between reference and Shimadzu Nexera X3

**Analyte:** Iohexol

**Measurand:** Iohexol in human serum

Characteristics	Value/Description
Measuring range	8.0 to 500 µg/ml
Limit of detection	<= 1.0 µg/ml
Limit of quantification	LLoQ <= 4 µg/ml with CV <= 15 % ULoQ >= 600 µg/ml with CV <= 15 %
Linearity	R >= 0.995 inside the measuring range No hook effect (evaluated up to 4,000 µg/ml)
Accuracy	5 % CV
Repeatability/Reproducibility	Inter-assay: <= 5 % (CI = 95 %) 20-day intra-assay: <= 3 % (CI = 95 %)

## Order information

Product No. (PN)	Product name
PN 001100	Nephrolyx IVDx Test Kit
PN 001200	Nephrolyx DNPx Digital Nephrology Basic Platform (1 year license)
PN 001310	Nephrolyx GFRx UHPLC Column
PN 001320	Nephrolyx GFRx UHPLC Guard Column

**Contact information:** [sales@nephrolyx.com](mailto:sales@nephrolyx.com)

## Contact

Nephrolyx GmbH  
Johann-Hittorf-Str. 8  
12489 Berlin  
Germany

[contact@nephrolyx.com](mailto:contact@nephrolyx.com)  
[www.nephrolyx.com](http://www.nephrolyx.com)



**For further information  
please contact:**

### **Dr Björn Neubacher**

Manager Sales & Business Development  
Phone +49 (0)174 1982066  
[sales@nephrolyx.com](mailto:sales@nephrolyx.com)

distributed in the US/Canada by:

**EAGLE BIOSCIENCES, INC.**

20A NW Blvd, Suite 112 Nashua, NH 03063  
Phone: 617-419-2019 FAX: 617-419-1110  
[www.EagleBio.com](http://www.EagleBio.com) • [info@eaglebio.com](mailto:info@eaglebio.com)



**EAGLE**  
BIOSCIENCES

**Nephrolyx**

May 2024