

### Liquichek Serum Indices

Monitor instrument response to detect pre-analytical HIL specimen interferences



### Improve Lab Performance

# Monitor the Ability of an Instrument to Detect Pre-Analytical Interferences

Hemolysis, Icterus and Lipemia (HIL) are the most commonly tested interferences that affect the integrity of patient samples in the clinical laboratory.<sup>1, 2</sup> Depending on the type and the degree of interferences, it may impact patient results when run with a colorimetric assay. Fortunately, most major chemistry platforms include an automated HIL detection feature that employs spectrophotometric detection to measure the level of common interferences. However, in lieu of automated detection, some laboratorians may visually inspect samples. In both scenarios, the level of interference is translated as indices (Hemolysis, Icterus and Lipemia index) according to Clinical and Laboratory Standards Institute (CLSI) Standard C56-A.

Liquichek Serum Indices is intended for use as part of laboratory practice to monitor the instrument response to detect Hemolyzed, Icteris or Lipemic (HIL) samples.

#### Similar to patient samples with human sourced material

Liquichek Serum Indices is prepared from human sourced material and is designed to monitor an instrument's response in detecting HIL specimens. This helps improve detection of pre-analytical errors and adds confidence to instrument performance.

#### Reduce human errors by eliminating manual inspections

Manual HIL interference detection involves subjective visual assessment from a trained laboratory technician that may result in longer turnaround times, errors or repeat testing. Monitoring the instrument response in HIL detection with Liquichek Serum Indices may help reduce hands-on time and ensure reliable results, enabling laboratory technicians to focus on other important tasks.

### Compare instrument responses among peer groups with the Unity Interlaboratory Program

An interlaboratory program can offer early awareness of shifts and trends to help avoid costly test repeats and unnecessary troubleshooting. Liquichek Serum Indices is integrated into the Unity Interlaboratory Program allowing access to instrument performance over time, peer group comparisons and easy-to-read reports.



#### **Regulators are taking note:** *ISO 15189:2012 includes expanded pre-examination procedures, and the IFCC\* is recommending standardized "Quality Indicators" in an effort to reduce the volume of pre-analytical errors.*<sup>1</sup>

Liquichek Serum Indices is a product that monitors instrument's response to detect HIL specimen interferences.

1. Green, S. F. The cost of poor blood specimen quality and errors in preanlytical processes. Clin BioChem. 2013;46:1175-9.

2. Salvagno, G., Lippi, G., Bassi, A., Poli, G., Guidi, G. Prevalence and type of pre-analytical problems for inpatient samples in coagulation laboratory. J Eval Clin Pract 2008;14:351–3.



### Liquichek Serum Indices

Liquichek Serum Indices monitors the instrument response to detect and flag HIL interference in a sample. Liquichek Serum Indices consists of three independent interferents (Hemolysis, Icterus, and Lipemia) and one non-interfered serum sample. The Hemolysis, Icterus, and Lipemia levels initiate instrument response based on increased hemoglobin, bilirubin, and lipoproteins respectively. The Non-interfered level is a human-based serum and is similar to a patient specimen. This product is unassayed, with target values and reference tables provided for convenience.<sup>3</sup> The reference tables include both quantitative indexes and qualitative interpretations compatible with major chemistry instrument platforms.

- Prepared from human source material
- 2 year shelf life at -20°C to -70°C
- 14 day open-vial stability at 2-8°C
- 14 day closed-vial stability at 2-8°C
- 28 day frozen aliquot at -20°C to -70°C

#### **Assay Sensitivity to HIL Interferences**

Interfering substances that remain undetected during pre-analytic inspection may impact the quality of results in the analytic testing cycle. The potentially negative consequences of inaccurate results must be prevented wherever possible. Some common assays with known sensitivity to HIL interferences are listed in the table below.

Hemolysis	Icterus	Lipemia
Amylase (AMY)	Amylase (AMY)	Aspartate Aminotransferase (AST)
Aspartate Aminotransferase (AST)	Creatinine (SC)	Direct Bilirubin (DBI)
Magnesium (MG)	Digoxin	Magnesium (MG)
Iron (FE)	Iron (FE)	Iron (FE)
Lactate Dehydrogenase (LDH)	γ-Glutamyl Transferase (GGT)	Lactate Dehydrogenase (LDH)
Direct Bilirubin (DBI)	Direct Bilirubin (DBI) Triglycerides (TG)	
Potassium (K)	Potassium/Sodium (K/Na)	Sodium (Na)

#### Indices

Hemolysis Icterus Lipemia Non-interfered

## Bio-Rad offers a comprehensive portfolio of products to cover all phases of clinical chemistry testing and data management



You can count on our consistent product quality to help assure reliability of patient results, along with expert technical support and customer educational programs.

#### **Ordering Information**

Cat #	Description	Cat #	Description
12012693	Hemolysis6 x 4 mL	12012697	Hemolysis MiniPak1 x 4 mL
12012694	Icterus6 x 4 mL	12012698	Icterus MiniPak1 x 4 mL
12012695	Lipemia6 x 4 mL	12012699	Lipemia MiniPak1 x 4 mL
12012696	Non-interfered6 x 4 mL	12012700	Non-interfered MiniPak1 x 4 mL

EQAS independent, external assessm Independent QC

Unity

An independent, external assessment of performance in comparison to your peers.

QCNet.com/eqas

Ongoing, proactive, unbiased daily QC that helps identify errors as they occur or begin to trend.

QCNet.com/independentqc

QC Data Management tools that help you create a strategy to reduce risk and streamline QC workflow.

QCNet.com/datamanagement



Bio-Rad Laboratories, Inc. For further information, please contact the Bio-Rad office nearest you or visit our website at www.qcnet.com/serumindices

Clinical Diagnostics Group Website www.bio-rad.com/diagnostics Australia +61 (2) 9914 2800 Austria +43 (0) 1 877 89 01 9 Belgium +32 (0) 3 710 53 00 Brazil +55 11 3065 7550 Canada +1 514 334 4372 China +86 21 6169 8500 Czech Republic +420 241 431 660 Denmark +45 44 52 100 Finland +358 9 804 22 00 France +33 (0) 1 47 95 60 00 Germany +49 (0) 89 31884 393 Greece +30 210 7774396 Hong Kong +85 2 2789 3300 Hungary +36 1 459 6190 India +91 124 4029300 Israel +972 03 963 6025 Italy +39 024 94 86 600 Japan +81 3 6361 7070 Korea +82 080 007 7373 Mexico +52 (55) 5488 7670 The Netherlands +31 (0)318 540 666 New Zealand +64 (9)415 2280 Norway +47 23 38 41 30 Poland +48 22 331 99 99 Portugal +351 21 47 27 700 Russia +7 495 721 1404 Singapore +65 6415 3170 South Africa +27 11 442 8508 Spain +34 91 490 6580 Sweden +46 844 98053 Switzerland +41 (0) 61 717 9555 Taiwan +886 (2) 2578-7189 Thailand (662) 651 8311 United Kingdom +44 (0)1923 471301

BIO-RAD and LIQUICHEK are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner.

© 2021 Bio-Rad Laboratories, Inc. Printed in USA QSD21-0053 Q-1652 Rev. 05/2021