

THE PERKINELMER HAND SANITIZER ANALYZER

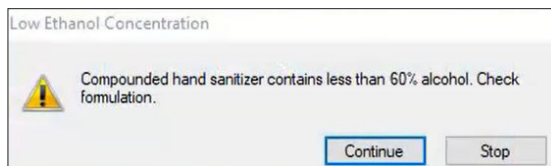


Recently the need for hygiene products have significantly increased, creating supply shortages. One of the most important of these is alcohol-based hand sanitizer. In order for alcohol-based hand sanitizers to truly decrease infectious agents on hands and other surfaces, they must meet specific defined criteria. It is known that hand sanitizers with insufficient alcohol concentration are not effective and could leave the user at higher risk of contracting and/or spreading infection. Furthermore the type of alcohol used is also critical to identify as some, such as Methanol/Wood alcohol, can be harmful to humans whereas ethanol or Isopropyl alcohol (IPA) are safer.

With many industries mobilizing to produce hand sanitizer for the first time, suppliers and distributors need quality control processes to ensure that the hand sanitizers produced or purchased are effective and safe.

The PerkinElmer Hand Sanitizer Analyzer is using an intuitive one-click operation and helps you:

- Determine if the product meets the requirement through a Pass/Fail information
- Maximize productivity with simple and rapid workflow driven verification



The PerkinElmer Hand Sanitizer Analyzer quickly and easily determines if the composition of the product meets guidelines for Hand Sanitizer released by WHO, USP or FDA.

Furthermore, with government and public concerns of counterfeit products, this analyzer can within seconds, verify efficacy of new products in the market.

The Analyzer is based on FT-IR technology and is self-installable, comes with easy step-by-step instructions and can be set up as easily as a printer. Remote support is available for any questions.



PerkinElmer Hand Sanitizer Analyzer.

For more information download the application brief at www.perkinelmer.com/handsanitizeranalyzer

You may also contact PerkinElmer directly via email: Asia Pacific: Sanitizer-APAC@Perkinelmer.com
Americas: Sanitizer-US@Perkinelmer.com, Europe, Middle East, Africa, India: Sanitizer-EMEA@Perkinelmer.com

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
P: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright © 2020, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

115630 (78724) PKI