

HAND SANITIZER TESTING SOLUTIONS



Label Your Brand With Confidence

From super simple alcohol screening, to accurate ingredients testing or multiple in-depth content analyses, varied hand sanitizer testing solutions are utilized depending on the needs of your business.

Find the PerkinElmer hand sanitizer testing solution that best fits your needs, while meeting FDA, WHO, USP and other guidelines or regulations. You will not only gain the right analyses for your operation; you will also gain support from the expert PerkinElmer team.



Several analytical testing solutions can provide you with accurate and precise results while optimizing your profits:

- Increased productivity with fast analysis
- Improved instrument uptime with our robust solutions
- Run samples immediately after installation with our turnkey analyzers and workflow solution guide

PerkinElmer Hand Sanitizer Analyzer – FT-IR	PerkinElmer Liquid Injection – GC/FID or GC with Confirmation	PerkinElmer Headspace Sampling – GC Method
Super Simple Screening	Accurate Label Claim	Multiple In-Depth Analyses
For end distributors, suppliers:	For alcohol producers, hand sanitizer manufacturers, suppliers, distributors QA/QC, and contract testing labs:	For hand sanitizer manufacturers, suppliers, distributors QA/QC, and contract testing labs:
<ul style="list-style-type: none"> • Simple affordable screening of ethanol and isopropanol content • Presence of harmful methanol down to 0.03 % • Fast, < 1 minute clear yes/ no screening of final product, determination if adhering to regulations • Difficulty level – low, a screening tool for anyone • No sample preparation required 	<ul style="list-style-type: none"> • Test raw materials (alcohol) for purity and contaminants per USP accurately and precisely • Identify and quantify impurities at detection limits defined by USP and FDA recently released criteria • Attain accurate Label Claim on your hand sanitizer • Avoid recalls or false positives using compound confirmation techniques 	<ul style="list-style-type: none"> • Expand your sample type with all the benefits of our Accurate Label Claim GC • Easy sample preparation for solid samples; know what's in your wipes • Avoid labor intensive and chemical laden extractions of solid samples for optimized productivity while reducing operating costs • Release your products faster to market

Please visit www.perkinelmer.com/handsanitizertesting

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.