



THE PERKINELMER CHLORINE ANALYZER

From keeping our drinking water safe to disinfecting our public transportation systems and hospitals, chlorine chemistry plays a critical role in the manufacturing of thousands of products we depend upon to remain healthy. Whether it's used in the manufacturing processes or in a final product, chlorine makes our world a safer place.

Due to the COVID-19 crisis, hygiene essentials are in such high demand, that there are now critical supply shortfalls for bleach and bleach-based wipe disinfectants. Underscoring the importance of disinfectants and sanitizer production, the U.S. Food and Drug Administration (FDA) released guidance for retail stores, restaurants, and food services during the pandemic to safeguard workers and consumers. Furthermore, the EPA's Emerging Pathogen Policy states that bleach products can combat COVID-19 when used as directed on hard, nonporous surfaces. Chlorine is the most important ingredient to produce bleach and bleach related products, however, due to chlorine's corrosive nature, manufacturers need analysis systems that are resistant to metal corrosion, while providing sensitivity and speed.

How does the PerkinElmer Chlorine Analyzer keep Chlorine manufacturers producing to meet the global need?

The **PerkinElmer Chlorine Analyzer** is a factory preconfigured and validated, turnkey gas chromatography (GC) solution, to obtain quick and reliable chlorine purity results, with a low cost of ownership.

- Exceed the requirements of the ASTM® E1746 "Standard Test Method for Sampling and Analysis of Liquid Chlorine for Gaseous Impurities"
- Run samples immediately upon installation
- Increase productivity with ultra-rugged, Chlorine corrosion resistant materials and venting capabilities for failure resistant operation
- Meet global demands by running 24/7

For more information visit: www.perkinelmer.com/chlorineanalyzer

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.