

MICROPLASTICS ANALYSIS WEBINAR SERIES

Optimizing Your Analytical Approach for Microplastics Identification

Each year, an estimated 9 million tons of plastic enters our lakes, rivers and oceans, eventually breaking down into the pollutants we now know as microplastics. With a lifespan of up to 450 years, microplastics materials can persist in the environment causing harm to marine life and entering the human food chain through seafood production.

This series of microplastics webinars is designed to help you overcome common testing challenges in your lab, understand the ecotoxicological effects of microplastics on marine life and recognize the importance of measures taken for quality assurance and control.

Who Would Benefit from Watching These Microplastics Analysis Webinars?

- Environmental Engineers
- Environmental Scientists
- Laboratory Managers and Directors

Quality Control Standards for Microplastic Analysis



Microplastics are recognized to be a global contaminant of concern, with the volume of both public attention and academic research on microplastics steadily increasing. This webinar details challenges faced in implementing contamination

controls, highlights the critical need for regulatory standards, and reviews the challenges in analyzing microplastics including extraction and quantification as well as the ability to compare results from different studies.

[WATCH NOW](#)

eXXpedition Round the World: Studying Plastic Pollution in our Global Ocean



In this webinar, the Science Lead of the all-women marine research group, eXXpedition, reviews the different scientific investigations and research operations carried out during their round-the-world expedition to study the flows of plastics from land into the environment and

analyze microplastics in surface water. The team used the PerkinElmer Spectrum Two™ FT-IR spectrometer which allowed them to analyze collected samples in near-real time with advanced FT-IR technology.

[WATCH NOW](#)

Emerging Pollutants and their Marine Ecotoxicological Impact



This webinar explores the ecotoxicological effects of microplastics and pharmaceuticals in marine ecosystems. These pollutants are a rising concern due to their wide-spread usage, limited removal by waste-water plants and frequent accumulation in tissues

of non-target species. Not only can microplastics induce slight cellular and physiological effects, they also interact in different ways with adsorbed chemical pollutants.

[WATCH NOW](#)

FT-IR as a 'Hub' Technique for Environmental Forensics and Engagement in Science



The application of FT-IR as a versatile analytical technique has increasingly played a central role in the investigative work in the Environmental sector, specifically in the Greenpeace Research Laboratories. This webinar provides an illustration of the ways in which FT-IR

instruments are used in laboratories, using examples of projects that have documented microplastics contamination, the complexity of marine microplastics as carriers of chemical burdens and the implications of the international trade in waste plastics for recycling and disposal.

[WATCH NOW](#)

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


PerkinElmer
For the Better

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

Copyright ©2021, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.