

A novel technique to qualify and quantify fine microplastics in surface water by coagulation and μ FTIR microscopy

○Yutaka Kameda

(Chiba Institute of Technology, Japan)

E-mail address: yutaka.kameda@it-chiba.ac.jp

Abstract

We propose an analytical method for fine micoplastics (MP) by using μ FTIR imaging, the Multivariate Curve Resolution (MCR) and the correlation analysis with their easy and quick sampling system. Coagulation was effective only for surface water samples with little suspended solids to remove matrices.

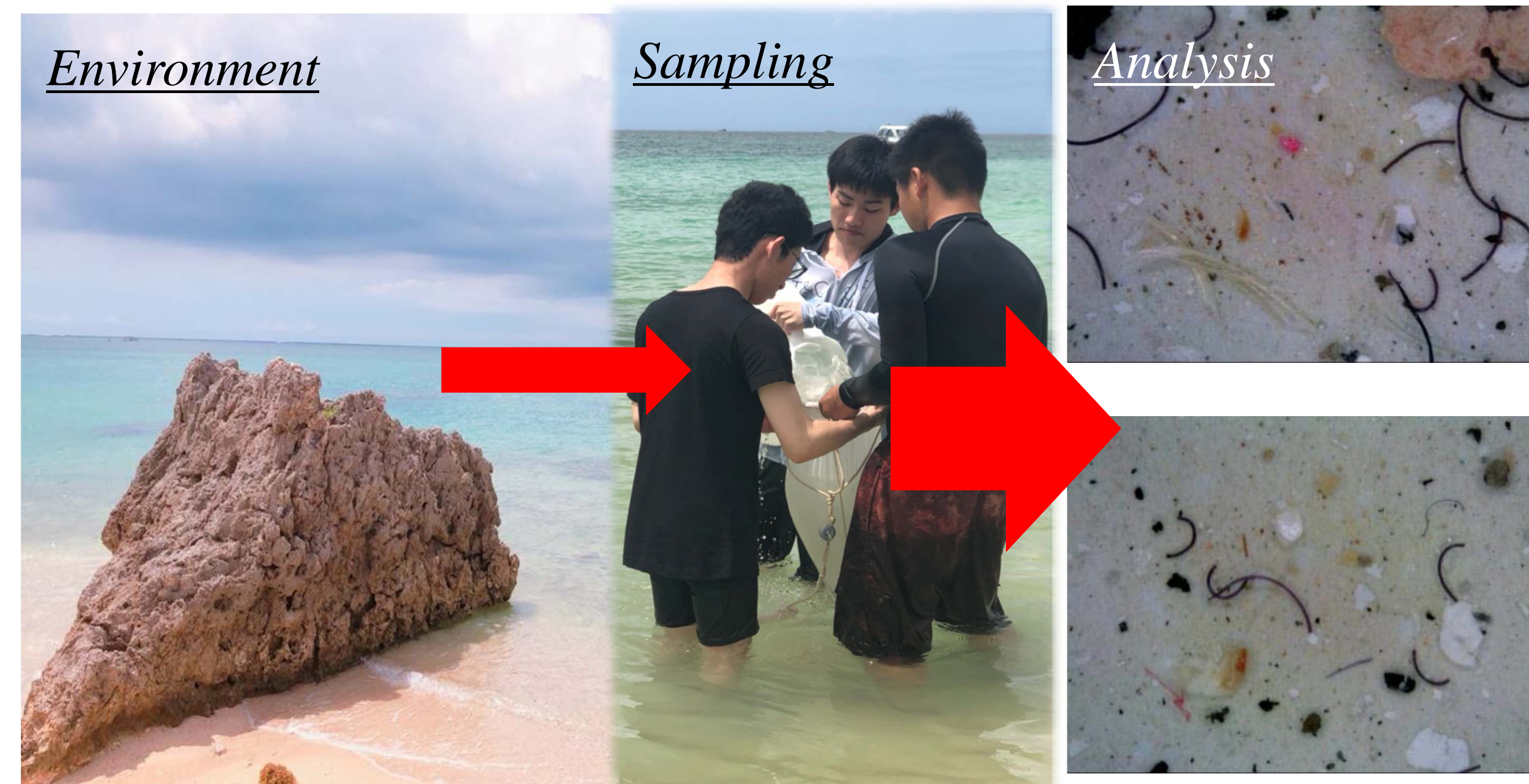


Fig. 1 Sampling and analysis of MP, especially fine MP, are difficult. Their identification by FTIR need traceability in their measurement.

Background

Fine MP (10 ~100 μ m) are growing concern with large MP (>100 μ m). The easy and quick sampling methods, the traceable analytical methods and the evaluation methods are needed.

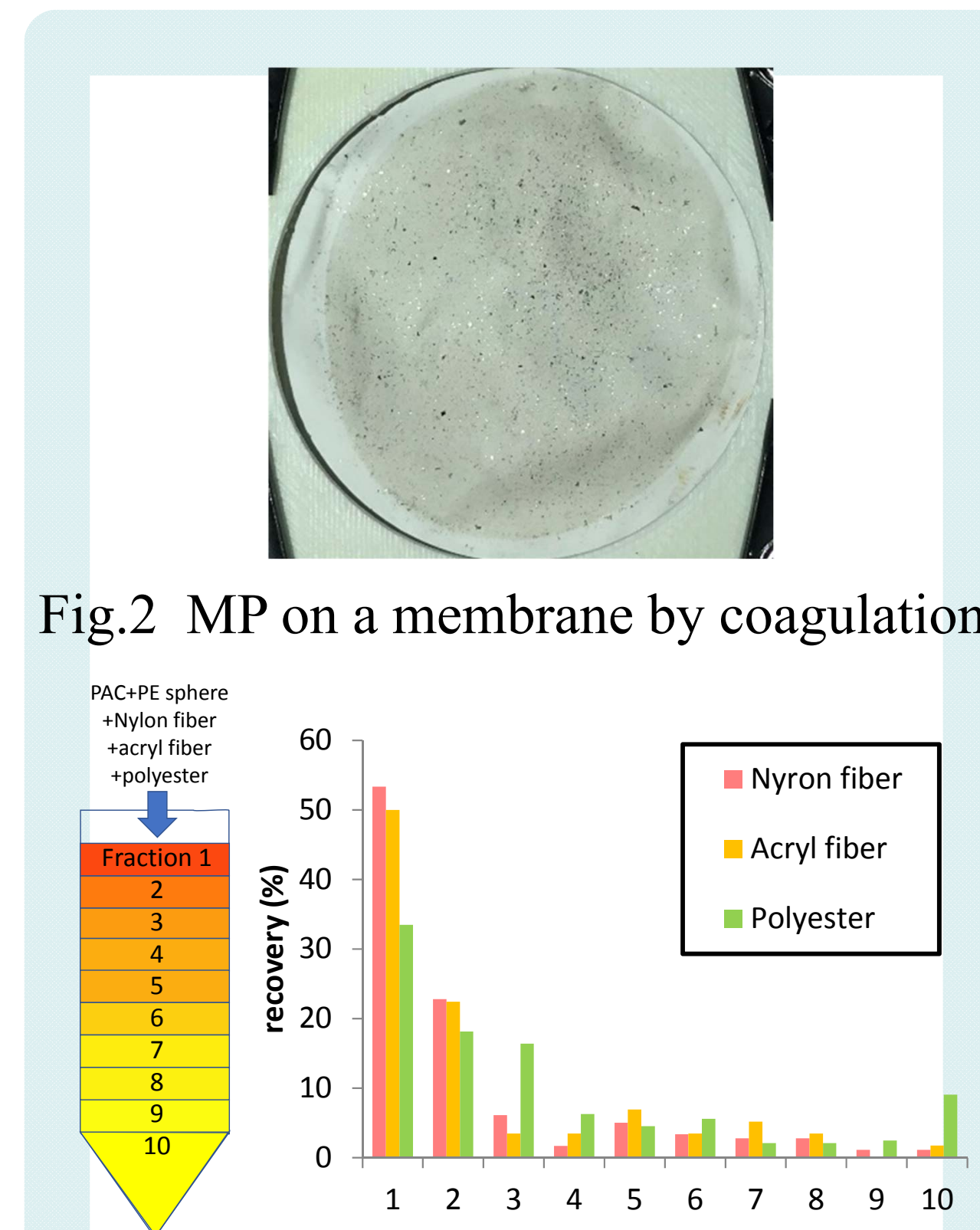


Fig.2 MP on a membrane by coagulation

Coagulation

We tried coagulation process after NaI density separation instead of dangerous H_2O_2 oxidation. Matrices such as algae, fine solid particles, plant debris could be removed easily and quickly (Fig. 2). Recovery tests (Fig. 3) using Milli Q with 30 μ m polyethylene particles and 100 μ m nylon, polyester and acryl fibers revealed good recovery rates. But cellulose fibers (toilet paper!) trapped the added MP and settled in raw wastewater.

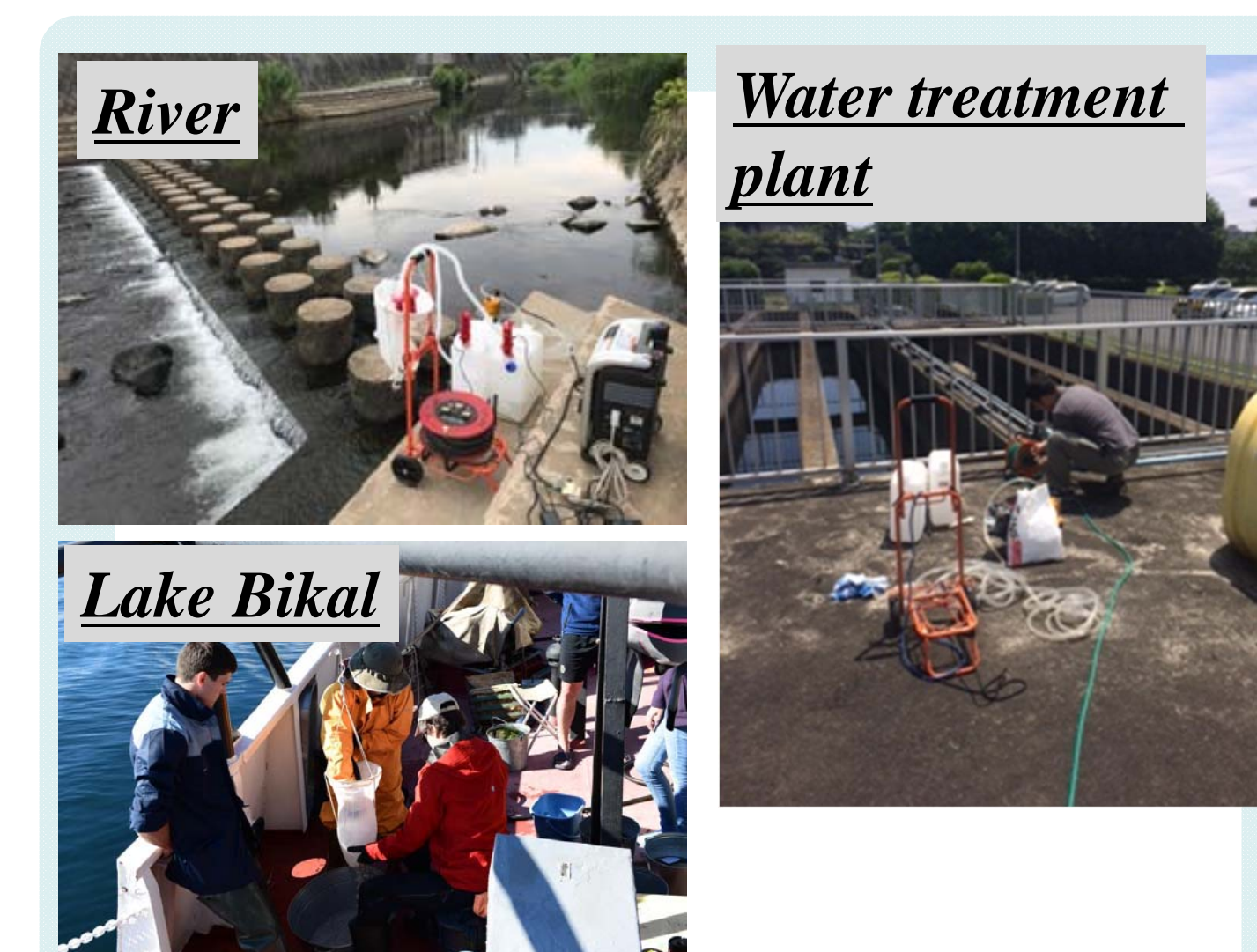


Fig. 4 The developed sampling system for 10 μ m MP are useful around the world.

Sampling System

Fig.4 shows a usage of a novel sampling system at various sites around the world. It takes 30 min to collect 1 m³ of water samples.

Automatic μ FTIR analysis

Fig. 5 shows an example of MP in surface water by automatic μ FTIR analysis. After sampling, MP were extracted on a PTFE membrane filter (2cm ϕ) after H_2O_2 oxidation and NaI density separation. MP images were conducted by μ FTIR (20% area/2hrs). Based on the spectrum data, MCR analysis can extract 10 major spectra. Minor spectrum can also be extracted by correlation analysis.

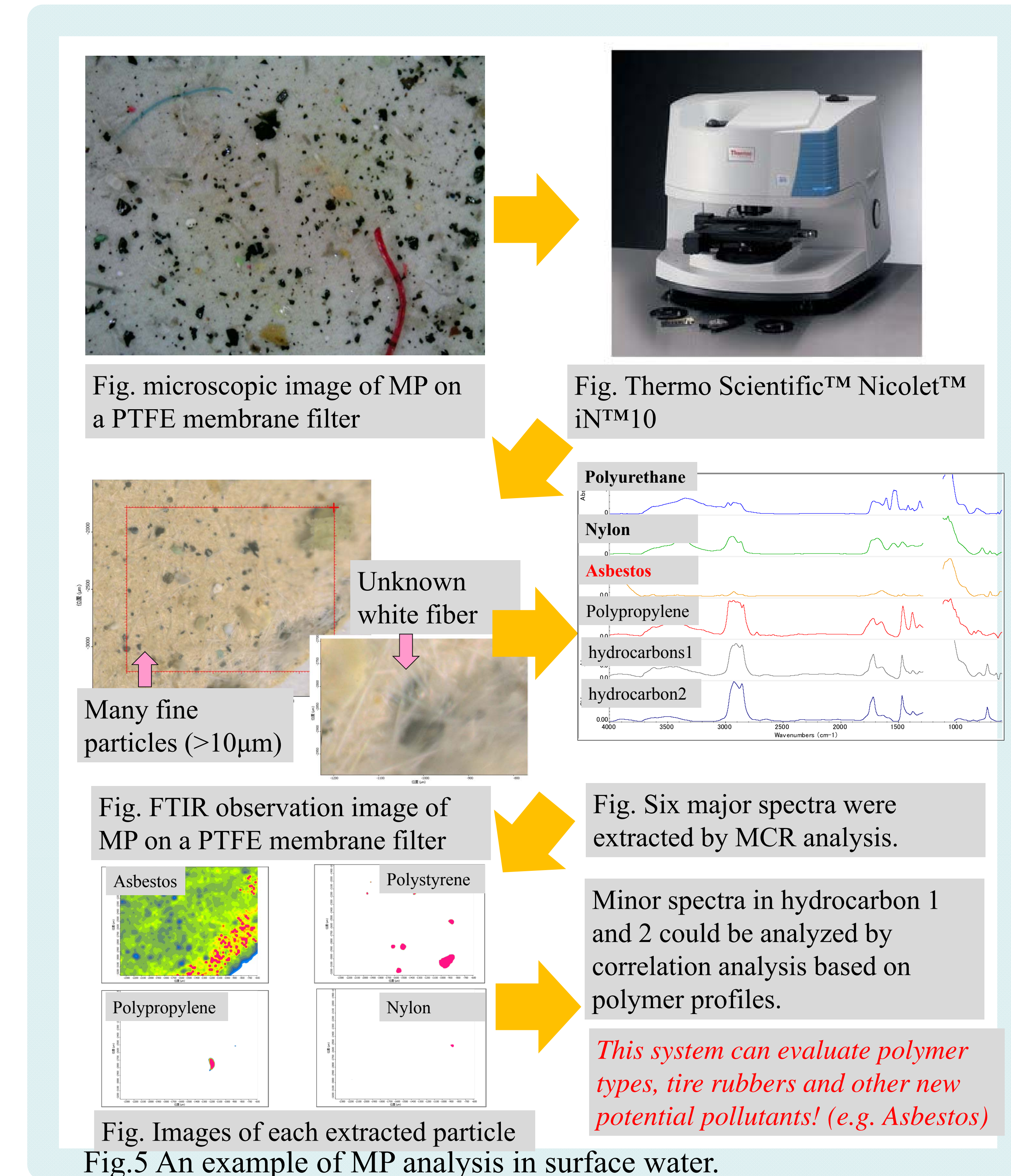


Fig.5 An example of MP analysis in surface water.