

An Unknown Screening Approach to Analyze Micro Pollutants Degradation by Disinfection Processes

Julia Quilitzk¹, Patricia Zerball-van Baar¹, Uwe Duennbier¹, Olaf Scheibner², Dipankar Ghosh³

¹Berliner Wasserbetriebe, Berlin, Germany

²Thermo Fisher Scientific, Bremen, Germany

³Thermo Fisher Scientific, San Jose, CA

ABSTRACT

In this study an unknown approach (suspect screening) in combination with a non-target screening was used to analyze micro pollutants degradation by disinfection processes. In addition 60 identified organic micro pollutants were quantified. The samples analyzed were collected at the wastewater treatment plant Ruhleben and include samples of treated wastewater (TW) before and after disinfection by UV radiation (UV) or performic acid (PA) treatment. The workflow from discovering a feature to the proposed structure using statistical tools is shown for one feature which was proposed to be triphenylphosphine oxide.

INTRODUCTION

Six wastewater treatment plants clean the wastewater from Berlin (Germany) and the surrounding regions. Treated wastewater flows into the rivers Spree and Havel. Ruhleben, the biggest wastewater treatment plant, is directly located in the city. For disinfection the last treatment step in Ruhleben is an UV radiation or performic acid treatment. These techniques can also be used in parallel if the waste water flow is too high. In August 2015 samples of the treated wastewater (TW) were taken before and after disinfection by UV radiation (UV) or performic acid (PA) treatment. This study shows a suitable way to identify "unknowns" or transformation products.

MATERIALS AND METHODS

Sample Preparation

Samples were directly injected onto the online-SPE system without dilution.

Liquid Chromatography

A 1 mL sample was injected onto the online-SPE system. For pre-concentration a C18 column, 2.1 × 20 mm with 12 μm particle size was used. For compound separation, a C18 column, 2.1 × 50 mm with 1.8 μm particle size was used. A gradient formed from water and methanol (B), both spiked with 0.1% formic acid, was ramped up from 2% B to 95% B in 6.7 minutes. The total chromatographic cycle time including online pre-concentration was 15 minutes.¹

Mass Spectrometry for Full Scan with variable Data Independent Acquisition

Mass spectrometric analysis was performed on a Thermo Scientific™ Q Exactive™ Focus hybrid quadrupole-Orbitrap™ mass spectrometer.

MS-parameter for screening and quantitative analysis in Full Scan Mode with variable Data Independent Acquisition (vDIA) or data dependent MS² (ddMS²):

for screening

Full scan (FS)

- ESI+ and ESI- separately
- R = 70,000; m/z 100–1000

variable Data Independent Acquisition (vDIA):

- R = 17,500
- CE = 20, 40, 60 eV
- vDIA segments: 5
- (m/z 100–205, 195–305, 295–405, 395–505, 495–1000)

data dependent MS² confirmation (ddMS²):

- R = 17,500
- CE = 20, 40, 60 eV
- Inclusion list with prioritized features after a statistical selection with Compound Discoverer

All resolution settings are provided in FWHM at m/z 200

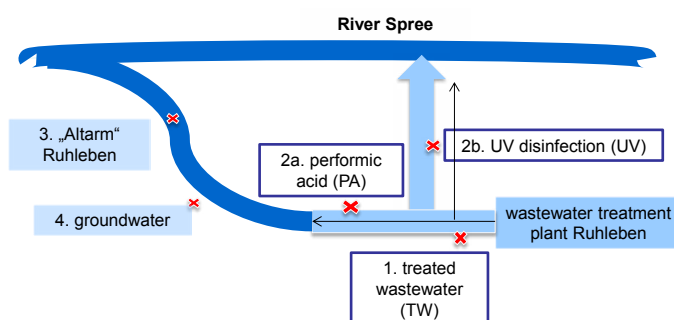
Note: vDIA method is not available in the United States of America

Data Analysis

Qualitative and quantitative analysis were performed with Thermo Scientific™ TraceFinder™ 3.3 software and Thermo Scientific™ Compound Discoverer™ 2.0 was used for the non-target screening approach.

SAMPLING

Figure 1. Schematic overview of the investigation area Berlin Ruhleben.



Investigation area

Figure 1 shows different sampling points in the area around the wastewater treatment plant Ruhleben. The following sampling points were in focus of the present study:

1. treated wastewater (TW) as control sample

The wastewater treatment plant has a throughput of around 247500 m³/day at dry weather. The wastewater treatment involves mechanical and biological purification steps. At first the water passes through a mechanically raked bar screen and six sand filters. This is followed by biological purification steps which include the removal of phosphate, nitrification and denitrification.²

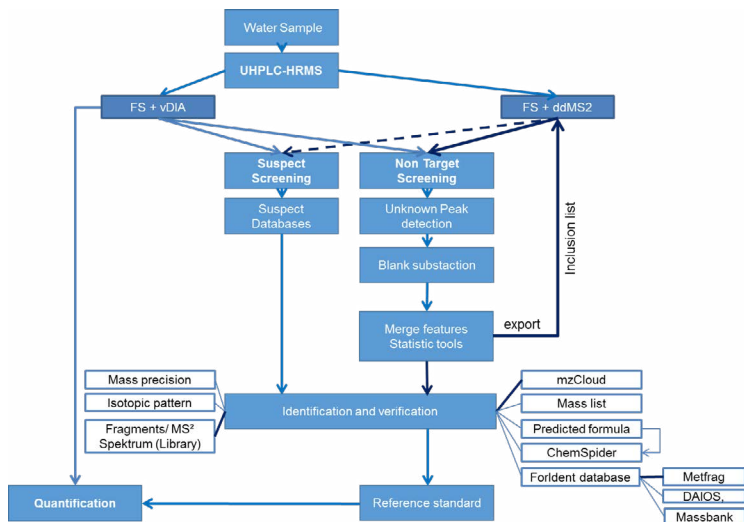
2. disinfection treatment steps (UV and PA) to evaluate any possible microbiological pollution degradation

2a. **Performic acid (PA):** For chemical disinfection performic acid is used in a concentration of 15–20ppm and the reaction time is about 8 minutes. The throughput amounts to 500 L/s.⁴

2b. **UV radiation (UV):** The UV radiation unit consists of a total of 648 radiation sources. The throughput is variable and ranges between 500 L/s and 1000 L/s.³

Workflow

Figure 2. Schematic overview of the workflow.



RESULTS - Suspect screening and quantification

Figure 2 shows a schematic diagram of the entire workflow carried out in this study. In the first step a suspect screening analysis was carried out with each sample (see fig.1). The database which was used contains a large number of compounds, e.g. pesticides, pharmaceuticals, industrial and household chemicals.

The suspect screening analysis resulted in a total of 381 suspects. By using confirmation criteria like mass precision and isotopic pattern match 192, 119 and 133 suspects could be confirmed in the control sample (TW) and the samples taken after UV or PA treatment, respectively. In addition 60 suspects could be quantified by reference standards. Table 1 shows the results of 17 selected compounds.

Table 1. Results of the suspect screening analysis and quantification.

Compound	Formula	used for/as	TW	UV	PA
			concentration [µg/L]		
Candesartan	C ₂₄ H ₂₀ N ₆ O ₃	treatment of hypertension	1.70	1.70 -	1.60 -
1H-Benzotriazole	C ₆ H ₅ N ₃	Corrosion inhibitor	3.90	4.00 -	3.40 ↓
2-Methylthiothiazole	C ₈ H ₇ NS ₂	vulcanization accelerator (metabolite)	0.47	0.55 -	x ↓
Bezafibrate	C ₁₉ H ₂₀ ClNO ₄	lipid-lowering agent	0.15	0.14 -	0.14 -
Bisoprolol	C ₁₈ H ₃₁ NO ₄	beta-blocker	0.43	0.51 -	0.25 ↓
Carbamazepine	C ₁₅ H ₁₂ N ₂ O	anticonvulsant	1.50	1.50 -	1.50 -
Diclofenac	C ₁₄ H ₁₁ Cl ₂ NO ₂	NSAID	1.40	0.86 ↓	1.1 ↓
10,11-Dihydro-10,11-dihydroxy carbamazepine	C ₁₅ H ₁₄ N ₂ O ₃	metabolite of carbamazepine	1.70	1.6 -	1.7 -
Carbendazim	C ₉ H ₉ N ₃ O ₂	fungicide	✓ []	✓ []	✓ []
Amisulpride	C ₁₇ H ₂₇ N ₃ O ₄ S	antipsychotic	✓ []	✓ []	✓ []
Irbesartan	C ₂₅ H ₂₈ N ₆ O	treatment of hypertension	✓ []	✓ []	✓ []
Clindamycin	C ₁₈ H ₃₃ ClN ₂ O ₅ S	antibiotic	✓ []	✓ []	x ↓
Sotalol	C ₁₂ H ₂₀ N ₂ O ₃ S	beta-blocker	0.18	0.17 -	0.09 ↓
Diethyltoluamide (DEET)	C ₁₂ H ₁₇ NO	insect repellents	0.21	0.28 -	0.12 ↓
Phenazone	C ₁₁ H ₁₂ N ₂ O	analgesic	✓ []	✓ []	✓ []
Oxazepam	C ₁₅ H ₁₁ ClN ₂ O ₂	benzodiazepine	✓ []	✓ []	✓ []
Primidone	C ₁₂ H ₁₄ N ₂ O ₂	anticonvulsant	0.29	0.32 -	0.35 -

✓ identified suspect but not quantified, x not detected

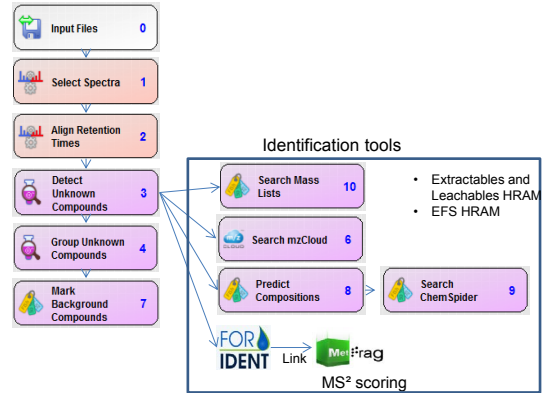
↓ concentration decrease relating to TW, - no concentration change relating to TW

The concentration of candesartan, carbamazepine, bezafibrate, 10,11-Dihydro-10,11-dihydroxy carbamazepine and primidone did not change after UV or PA disinfection. Whereas the concentration of 1H-benzotriazole, 2-methylthiothiazole, bisoprolol, sotalol, diclofenac and diethyltoluamide decreased after treatment with PA. In addition the Diclofenac concentration also decreased significantly after UV treatment. That is well-known for diclofenac.

RESULTS - Non-Target Screening

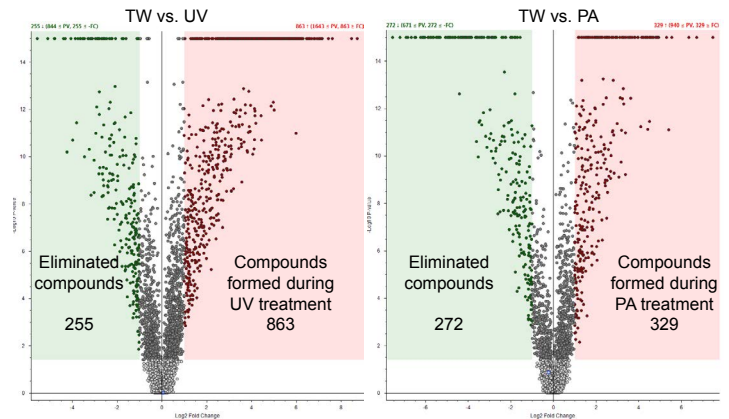
To identify further unknown compounds an unknown approach with Compound Discoverer 2.0 was conducted.

Figure 3. Unknown approach with Compound Discoverer 2.0 (workflow).



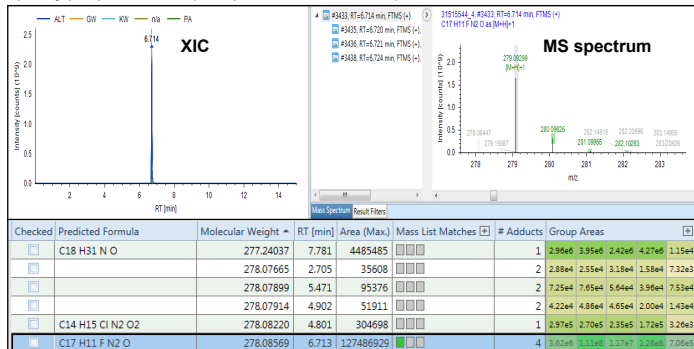
Compound Discoverer 2.0 (figure 3) is used to detect unknown compounds via elemental composition predictions. The software automatically subtracts the background (blanks), performs mzCloud and ChemSpider library searchings and scores the compounds based on defined patterns. In addition the database ForIdent which includes only water relevant data was used in this approach.

Figure 4. Unknown approach with Compound Discoverer 2.0 (workflow).



Statistic tools are useful to identify interesting features. In figure 4 the compounds formed after treatment with UV or PA are shown in red. At first glance it seems that more „new“ features were formed after UV treatment. A total of 15154 compounds were detected in all samples. After filtering the data by UV/TW >10 or PA/TW >10 the number of unique features decreased to 863 and 329 in the UV and PA samples, respectively. Only those features were in focus of this study.

FIGURE 5. Screenshot of compound m/z 278.08569 with masslist search hit triphenylphosphine oxide (Compound Discoverer 2.0).



Identification results

Checked	ΔMass [Da]	ΔMass [ppm]	Formula	Molecular Weight	RT [min]	Structure	Name
<input checked="" type="checkbox"/>	0.00036	1.31	C18 H15 O P	278.08605	0.000		Triphenylphosphine oxide

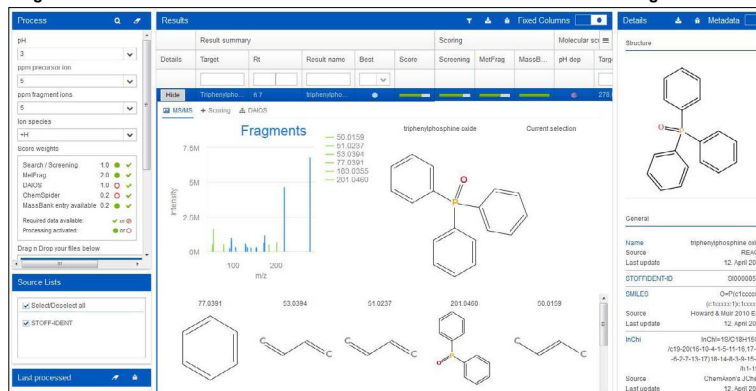
Checked	ΔMass [Da]	ΔMass [ppm]	Scan Number	mcCloud ID	Formula	Structure	Molecular Weight	Match	Best Match	Name
<input checked="" type="checkbox"/>	0.00027	0.98	1487	92	C18 H15 O P		278.08605	71.5	71.5	Triphenylphosphine oxide

Figure 5 shows a selected compound with the monoisotopic mass 278.08569. Besides [M+H]⁺ the adducts [M+Na]⁺, [2M+H]⁺ and [2M+Na]⁺ were also detected. This compound was only found in the UV sample. The masslist search resulted in triphenylphosphine oxide (TPPO) which is a waste product in various chemical reactions e.g. Wittig reaction. A library score of 71.5% was given by mzcloud.

FOR-IDENT is an online database especially for compounds relevant in water analysis. 9500 compounds are currently registered and the database is continuously updated and extended. A feature list with approximately 5000 features including their exact masses and retention times can be uploaded. The main criteria used for compound identification include exact masses (max. tolerance set at Δ 5 ppm), fragment ions (Δ 5 ppm), the ion species (+H) and the retention time index (RTI). If MS2-data are available the linked in MetFrag tool can be used additionally to further prioritize suggested compounds via *in silico* fragmentation.

In case of the compound with m/z 279.09349 the structure with the highest score is triphenylphosphine oxide (see figure 6). For final confirmation a reference standard needs to be checked.

Figure 6. Forident-Hit for the mass 279.09349 with ddMS2- information from MetFrag



CONCLUSION

The present study is a combination of quantification and screening. With the suspect screening a high number of suspect could be identified and verified by reference standards. Whereby the behaviour of the different compounds could be observed. The quantified compounds showed no significant decrease of concentration by the disinfection step UV with the exception of diclofenac. With PA treatment often the concentration decreases. With statistical tools in the non-target approach the data set can be reduced significantly to figure out relevant features. For identification beside isotopic pattern and formula prediction a pattern with databases and libraries are feasible. The procedure of the non-target workflow is shown from Pkfinding to identifying with an example. To identify „new“ unknowns results it is a long and difficult way to discover.

REFERENCES

1. Wode et al. JCA **2012**, 1270, 118–126.
2. <http://www.bwb.de/content/language1/html/981.php>, access date: May 2016
3. Regina Gnirss; Ulrike Franzke, http://kompetenz-wasser.de/fileadmin/user_upload/pdf/veranstaltungen/OXERAM_Info-Tag/20110525_Infotag_Oxeram_BWB.pdf, 2011, access date: May 2016
4. Veronika Gronau, Optimierung und Sicherung der Datenqualität von STOFF-IDENT; Wasser 2016; Jahrestagung der wasserchemischen Gesellschaft, 2.5.-4.5.2016, Bamberg, 449-453

ACKNOWLEDGEMENTS

The authors thank the BMBF for further funding in the continuous strategy development via the project initiative „FOR-IDENT“ (support code: 02WRS1354B).



www.thermofisher.com

©2016 Thermo Fisher Scientific Inc. All rights reserved. ISO is a trademark of the International Standards Organization. All other trademarks are the property of Thermo Fisher Scientific and its subsidiaries. This information is presented as an example of the capabilities of Thermo Fisher Scientific products. It is not intended to encourage use of these products in any manners that might infringe the intellectual property rights of others. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.

Africa +43 1 333 50 34 0
Australia +61 3 9757 4300
Austria +43 810 282 206
Belgium +32 53 73 42 41
Brazil +55 11 2730 3006
Canada +1 800 530 8447
China 800 810 5118 (free call domestic)
 400 650 5118

Denmark +45 70 23 62 60
Europe-Other +43 1 333 50 34 0
Finland +358 10 3292 200
France +33 1 60 92 48 00
Germany +49 6103 408 1014
India +91 22 6742 9494
Italy +39 02 950 591

Japan +81 6 6885 1213
Korea +82 2 3420 8600
Latin America +1 561 688 8700
Middle East +43 1 333 50 34 0
Netherlands +31 76 579 55 55
New Zealand +64 9 980 6700
Norway +46 8 556 468 00

Russia/CIS +43 1 333 50 34 0
Singapore +65 6289 1190
Sweden +46 8 556 468 00
Switzerland +41 61 716 77 00
Taiwan +886 2 8751 6655
UK/Ireland +44 1442 233555
USA +1 800 532 4752
 PN64758-EN 0616S

