

使用超高灵敏度三重四极杆质谱仪实现水中药物和个人护理用品 (PPCP) 的高通量检测

作者

杜伟、李建中、郭启雷、薄涛
安捷伦科技（中国）有限公司

摘要

本文采用安捷伦三重四极杆液质联用系统，建立了测定环境水中药物和个人护理用品 (PPCP) 的分析方法。文中利用直接进样方式，应用动态 MRM 采集模式实现了同时对环境水中 377 种 PPCP 快速准确的筛查和定量分析，多数化合物的检测限 (LOD) 低于 10 ng/L，无需进行样品富集。该方法线性关系良好，在添加浓度为 100 ng/L 时，超过 63% 的化合物的回收率在 80%-120% 之间，超过 78% 的化合物的 RSD < 20%，重复性和回收率良好，完全满足环境领域监测需求。

前言

药物和个人护理用品 (Pharmaceutical and Personal Care Products, PPCP) 作为一类新兴污染物日益受到人们的关注。PPCP 种类繁多, 包括各类抗生素、止痛药、降压药、避孕药、催眠药、减肥药、发胶、染发剂、香水和杀菌剂等, 人们的使用是 PPCP 进入环境的主要方式, 它通常以痕量浓度存在于地表水中, 对环境、野生动植物和人类的健康带来负面影响。

环境水中的 PPCP 浓度很低, 通常处于 ng/L 水平。采用常规的三重四极杆液相色谱与质谱联用系统进行检测时, 通常需要通过固相萃取 (SPE) 对样品进行富集。SPE 需要大量样品, 操作繁琐、耗时长, 成为限制检测通量和准确性的因素。随着超高灵敏度 Agilent 6495 三重四极杆质谱仪结合安捷伦喷射流离子源的出现, 可以采用直接进样的方式检测环境水中低至 ng/L 浓度水平的 PPCP, 简化了前处理过程, 显著提高了检测通量。

本文以直接进样方式, 采用 Agilent 1290 Infinity II 液相色谱/6495 三重四极杆液质联用系统, 建立了同时检测 377 种 PPCP 类化合物的可靠定量方法。包括抗生素类 104 种、兴奋剂类 29 种、激素类 46 种、精神类 48 种、解热镇痛类 28 种、抗过敏类 13 种、心血管类 20 种、农药 48 种以及其它类 41 种。该方法具有高通量、高灵敏度和覆盖范围广等特点, 适合于环境行业 PPCP 高通量的筛查和检测。

实验部分

试剂和样品

所有试剂均为 HPLC-MS 级。乙腈购自 Honeywell。超纯水产自 Milli-Q 纯水系统 (Merck KGaA)。5 mol/L 乙酸铵溶液购自 Fluka (09691-250ML)。乙酸购自 Aldrich (338828-25ML)。PPCP 标准品由外部合作方提供。实际水样取自于北京某湖泊及其周边水域。

仪器和设备

实验采用 Agilent 1290 Infinity II 液相色谱进行分离, 其由下列部件组成:

- Agilent 1290 Infinity II 二元高速泵 (G7120A)
- Agilent 1290 Infinity II 高通量进样器 (G7167B)
- Agilent 1290 Infinity II 多柱柱温箱 (G7116B)

采用配有安捷伦喷射流 (AJS) 电喷雾离子源和 iFunnel 离子传输聚焦技术的 Agilent 6495 三重四极杆液质联用系统, 在正离子模式下进行检测。

采用 Agilent MassHunter 工作站软件 (B.07 版本) 进行数据采集和分析。

标准溶液配置

PPCP 标准品储备液以乙腈配制, 各化合物的浓度均为 10 mg/L。以 Milli-Q 纯水配制各化合物的混合标准溶液, 校准曲线浓度范围为 10-10000 ng/L。浓度为 100 ng/L 的加标样品由实际水样中添加混合标准溶液配制。

样品前处理

实际水样采用 0.22 μ m 滤膜过滤后直接进样 20 μ L 进行分析检测。

表 1. 用于正离子模式方法的液相色谱条件

色谱柱	ZORBAX Eclipse Plus C18, 2.1 × 100 mm, 1.8 μm (部件号: 959758-902)	
柱温	40 °C	
进样量	20 μL	
流动相	A) 5 mM 乙酸铵 + 0.02% 乙酸的水溶液 B) 乙腈	
流速	0.3 mL/min	
梯度程序	时间 (min)	B(%)
	0	5
	0.5	5
	11	100
	15	100
停止时间	15 min	
后运行时间	3 min	

表 2. 用于负离子模式方法的液相色谱条件

色谱柱	ZORBAX Eclipse Plus C18, 2.1 × 100 mm, 1.8 μm (部件号: 959758-902)	
柱温	40 °C	
进样量	20 μL	
流动相	A) 0.005% 乙酸水溶液 B) 乙腈	
流速	0.3 mL/min	
梯度程序	时间 (min)	B(%)
	0	5
	0.5	5
	6	100
	9	100
停止时间	15 min	
后运行时间	3 min	

表 3. 质谱离子源条件

离子模式	正离子模式/负离子模式
干燥气温度	250 °C (正) /200 °C (负)
干燥气流速	16 L/min (正) /12 L/min (负)
鞘气温度	380 °C
鞘气流速	12 L/min
雾化器压力	40 psi
毛细管电压	3000 V
喷嘴电压	0 V (正) /2000 V (负)
MS ₁ 和 MS ₂ 分辨率	单位分辨率
MRM 参数	如表 4 和表 5 所示

表 4. 正离子模式下化合物 MRM 离子对参数列表

化合物名称	母离子 (m/z)	子离子 (m/z)	碰撞能量 (V)	保留时间 (min)	化合物名称	母离子 (m/z)	子离子 (m/z)	碰撞能量 (V)	保留时间 (min)
10,11-dihydro-10-hydroxycarbamazepine	255.1	237	20	4.51	Lidocaine	235.1	86.2	8	4.03

化合物名称	母离子 (m/z)	子离子 (m/z)	碰撞能量 (V)	保留时间 (min)	化合物名称	母离子 (m/z)	子离子 (m/z)	碰撞能量 (V)	保留时间 (min)
10,11-dihydro-10-hydroxycarbamazepine	255.1	194.1	4	4.51	Lidocaine	235.1	58.1	32	4.03
2-Aminobenzimidazole	134.1	92	25	2.60	Lincomycin	407.2	359	15	3.23
2-Aminobenzimidazole	134.1	65.1	35	2.60	Lincomycin	407.2	126	30	3.23
2-aminoflubendazole	256.1	123	40	4.84	LMG	331.2	239.1	30	10.90
2-aminoflubendazole	256.1	95	40	4.84	LMG	331.2	223.1	55	10.90
2-Mercaptobenzimidazole	151	93	15	3.63	Loratadine	383.2	337.1	20	8.96
2-Mercaptobenzimidazole	151	65	20	3.63	Loratadine	383.2	267.1	40	8.96
2-Methyl-4(5)-nitroimidazole	128.1	82	16	2.04	Lorazepam	321	275.1	12	6.11
2-Methyl-4(5)-nitroimidazole	128.1	42	36	2.04	Lorazepam	321	229.2	28	6.11
2-Methyl-5-nitroimidazole	128.1	82	16	3.50	Maduramycin	934.4	647.3	16	12.38
2-Methyl-5-nitroimidazole	128.1	42	36	3.50	Maduramycin	934.4	629.4	20	12.38
2-Thiouracil	129	112	12	1.19	Malathion	331	127	5	8.39
2-Thiouracil	129	70	17	1.19	Malathion	331	99	10	8.39
3-Hydroxycarbofuran	238	181	5	4.37	MDA	180.1	163.2	4	3.32
3-Hydroxycarbofuran	238	163	10	4.37	MDA	180.1	105.2	20	3.32
4-Aminoantipyrine	204.1	159	8	4.05	MDEA	208.1	163.1	4	3.71
4-Aminoantipyrine	204.1	83	12	4.05	MDEA	208.1	77.1	44	3.71
4-Isopropylaminoantipyrine	246.2	153.2	8	5.02	MDMA	194.1	163.1	8	3.46
4-Isopropylaminoantipyrine	246.2	56.2	20	5.02	MDMA	194.1	77.1	40	3.46
4-Nitroimidazole	114.1	97	15	1.47	Mebendazole	296.1	264.1	18	5.98
4-Nitroimidazole	114.1	68	25	1.47	Mebendazole	296.1	105	35	5.98
5-Chloro-1-methyl-4-nitroimidazole	162.1	145	16	3.56	Mefenamic acid	242.1	224	16	7.59
5-Chloro-1-methyl-4-nitroimidazole	162.1	116	18	3.56	Mefenamic acid	242.1	208	36	7.59
5-Hydroxymebendazole/5-Hydroxyl-mebendazole	298.2	266.1	19	5.20	Megestrol acetate	385.2	325.2	8	8.70
5-Hydroxymebendazole/5-Hydroxyl-mebendazole	298.2	160	36	5.20	Megestrol acetate	385.2	267.2	12	8.70
5-Hydroxy-Thiabendazole	218	176	17	5.93	Melengestrol acetate	397.2	337.3	8	8.88
5-Hydroxy-Thiabendazole	218	148	25	5.93	Melengestrol acetate	397.2	279.2	20	8.88
5-Nitrobenzimidazole	164.1	118	22	4.10	Meperidine	248	220.2	16	4.49
5-Nitrobenzimidazole	164.1	91	40	4.10	Meperidine	248	174.1	16	4.49
6-Acetylmorphine	328.1	211.1	24	3.31	Meprobamate	219	158.1	0	4.65
6-Acetylmorphine	328.1	165.1	44	3.31	Meprobamate	219	97	12	4.65
Acebutolol	337.2	116	16	3.92	Mercaptodimethur	226.3	169	7	7.53
Acebutolol	337.2	56	40	3.92	Mercaptodimethur	226.3	121	20	7.53
Acetaminophen	152	110	12	2.54	Metaproterenol/Orciprenaline	212.1	194.1	6	1.54
Acetaminophen	152	65.1	32	2.54	Metaproterenol/Orciprenaline	212.1	152	14	1.54
Alanycarb	422	199	15	8.40	Metformin	130.1	71.1	24	0.62
Alanycarb	422	142	20	8.40	Metformin	130.1	60	12	0.62
Albendazole	266.1	234.1	16	6.69	Methadone	310.2	265.2	8	6.19
Albendazole	266.1	191	34	6.69	Methadone	310.2	104.9	28	6.19

化合物名称	母离子 (m/z)	子离子 (m/z)	碰撞能量 (V)	保留时间 (min)	化合物名称	母离子 (m/z)	子离子 (m/z)	碰撞能量 (V)	保留时间 (min)
Albendazole-2-aminosulfoxide	240.1	198.1	16	3.55	Methamphetamine	150.1	119.1	8	3.39
Albendazole-2-aminosulfoxide	240.1	133	30	3.55	Methimazole/Tapazole/TAP	115	88	16	1.53
Albendazole-Sulfone	298.1	266.1	16	4.82	Methimazole/Tapazole/TAP	115	57	19	1.53
Albendazole-Sulfone	298.1	159	39	4.82	Methiocarb	226	169	5	7.42
Albendazole-Sulfoxide	282.1	240.1	7	4.23	Methiocarb	226	121	10	7.42
Albendazole-Sulfoxide	282.1	208	21	4.23	Methiocarb sulfone	258.2	201	13	5.01
Albuterol	240.1	222.2	0	2.39	Methiocarb sulfone	258.2	121.7	15	5.01
Albuterol	240.1	148	12	2.39	Methiocarb sulfoxide	242.2	185	17	4.11
Alclomethasone dipropionate	521.2	301	10	8.53	Methiocarb sulfoxide	242.2	122	17	4.11
Alclomethasone dipropionate	521.2	279	10	8.53	Methomyl	163	106	5	3.50
Aldicarb	116	89	5	5.37	Methomyl	163	88	5	3.50
Aldicarb	116	70	5	5.37	Methomyl-oxime	106	88	7	3.48
Aldicarb sulfone	223	148	5	3.26	Methomyl-oxime	106	58	10	3.48
Aldicarb sulfone	223	86	15	3.26	Methotrexate	455.1	308.2	16	2.88
Aldicarb sulfoxide	207	132	5	2.74	Methotrexate	455.1	175.1	36	2.88
Aldicarb sulfoxide	207	89	5	2.74	Methylphenidate	234.1	84.2	20	4.17
Aminonide	503.2	338.9	10	8.54	Methylphenidate	234.1	56.2	52	4.17
Aminonide	503.2	321	14	8.54	Methylprednisolone	375.2	357.1	6	5.83
Aminocarb	209.1	152	10	5.57	Methylprednisolone	375.2	161.1	20	5.83
Aminocarb	209.1	137	20	5.57	Methylprednisolone 21-acetate	417.2	399.2	6	6.99
Amino-Mebendazole/ Mebendazole-amine(HMEB)	238	105	25	4.58	Methylprednisolone 21-acetate	417.2	253.2	18	6.99
Amino-Mebendazole/ Mebendazole-amine(HMEB)	238	77	40	4.58	Methylthiouracil/6-Methyl-2-thiouracil	143	126	14	1.59
Aminophenazone	232	113	10	4.54	Methylthiouracil/6-Methyl-2-thiouracil	143	84	15	1.59
Aminophenazone	232	111	10	4.54	Metolcarb	166.1	109	5	2.73
Amitriptyline	278.2	202.2	68	6.11	Metolcarb	166.1	94	35	2.73
Amitriptyline	278.2	91	24	6.11	Metoprolol	268.2	76.9	56	4.05
Amitriptyline metabolite	294.2	276.2	8	4.56	Metoprolol	268.2	56.2	24	4.05
Amitriptyline metabolite	294.2	214.9	48	4.56	Metronidazole	172.1	128	12	2.82
Amphetamine	136.1	119.1	4	3.22	Metronidazole	172.1	82	26	2.82
Amphetamine	136.1	91.1	12	3.22	Mevastatin	391.2	185.1	8	8.86
Aripiprazole	448.1	285.1	24	6.75	Mevastatin	391.2	159.1	24	8.86
Aripiprazole	448.1	98.2	40	6.75	MG	329.2	314.2	38	7.18
Astemizole	459.3	218.2	20	6.55	MG	329.2	285.1	42	7.18
Astemizole	459.3	135.1	40	6.55	m-Hydroxybenzoylecgonine	306.1	168.1	12	3.32
Atenolol	267.2	145.1	24	2.50	m-Hydroxybenzoylecgonine	306.1	65.2	72	3.32
Atenolol	267.2	74	20	2.50	Miconazole	415	159	40	9.97
Atorvastatin	559.2	440.2	20	6.95	Miconazole	415	69	20	9.97

化合物名称	母离子 (m/z)	子离子 (m/z)	碰撞能量 (V)	保留时间 (min)	化合物名称	母离子 (m/z)	子离子 (m/z)	碰撞能量 (V)	保留时间 (min)
Atorvastatin	559.2	250.2	40	6.95	MNZOH/Hydroxy metronidazole/2-hydroxy-methyl-5-nitroimidazol	188.1	126	15	2.21
Atrazine	216.1	174.1	8	6.51	MNZOH/Hydroxy metronidazole/2-hydroxy-methyl-5-nitroimidazol	188.1	123	10	2.21
Atrazine	216.1	68.2	36	6.51	Modafinil	296.1	129.2	8	5.17
Avermectin	895.5	751.4	45	11.12	Mometasone furoate	521.1	503	4	8.66
Avermectin	895.5	449.2	50	11.12	Mometasone furoate	521.1	263	24	8.66
Azaperol	330	149	25	4.73	Monensin-NH4	688.5	635.4	12	12.65
Azaperol	330	121	30	4.73	Monensin-NH4	688.5	461.3	23	12.65
Azaperone	328	165	7	5.30	Monoethylglycinexylidide	207.1	122.2	8	3.37
Azaperone	328	121	21	5.30	Monoethylglycinexylidide	207.1	58.2	4	3.37
Azlocillin	462	246	10	3.53	Montelukast	586.2	422.1	20	10.32
Azlocillin	462	218	20	3.53	Montelukast	586.2	278.1	28	10.32
Beclomethasone	409.2	391.1	6	6.13	Morphine	286.1	152.2	64	2.02
Beclomethasone	409.2	146.9	30	6.13	Morphine	286.1	128	68	2.02
Beclomethasone dipropionate	521.2	503	4	9.06	Naftifine	288.2	141.1	20	8.61
Beclomethasone dipropionate	521.2	319	10	9.06	Naftifine	288.2	117.1	20	8.61
Bendiocarb	224	167	5	6.14	Nalidixic Acid	233	215	10	5.91
Bendiocarb	224	109	10	6.14	Nalidixic Acid	233	187	20	5.91
Benfuracarb	411	252	10	9.90	Nandrolone/Nortestosterone	275.2	239.1	13	6.67
Benfuracarb	411	195	20	9.90	Nandrolone/Nortestosterone	275.2	109	24	6.67
Bentazon	241	184	8	6.57	Naproxen	231	185	10	6.43
Bentazon	241	57	20	6.57	Naproxen	231	141	40	6.43
Benzimidazole	119.1	92	30	2.65	Narasin	787.6	531.3	52	13.08
Benzimidazole	119.1	65.1	35	2.65	Narasin	787.6	431.2	56	13.08
Benzoylcegonine	290.1	168.2	16	3.55	Nifedipine	347.1	315.2	0	7.01
Benzoylcegonine	290.1	77	64	3.55	Nifedipine	347.1	195.1	36	7.01
Betamethasone valerate	477.3	354.9	4	8.14	Nifedipine oxidized	345.1	284	24	6.92
Betamethasone valerate	477.3	278.8	14	8.14	Nifedipine oxidized	345.1	268.1	24	6.92
Betamethasone dipropionate	505.2	318.9	10	8.79	Nigericin M+NH4	742.5	675.3	25	13.82
Betamethasone dipropionate	505.2	278.9	12	8.79	Nigericin M+NH4	742.5	657.2	25	13.82
Bifonazole	311.2	243.1	10	8.66	Nimorazole	227.1	114.1	10	3.56
Bifonazole	311.2	165.1	40	8.66	Nimorazole	227.1	100.1	20	3.56
Brompheniramine	319.1	274	20	5.20	Norfentanyl	233.1	84	12	3.74
Brompheniramine	319.1	167.1	40	5.20	Norfentanyl	233.1	55	40	3.74
Budesonide	431.2	413.1	6	7.26	Norfluoxetine	296.1	134.1	0	6.00
Budesonide	431.2	146.9	30	7.26	Norfluoxetine	296.1	30	8	6.00
Buprenorphine	468.3	84.2	48	7.51	Normeperidine	234.1	160.3	8	4.41
Buprenorphine	468.3	55.1	52	7.51	Normeperidine	234.1	91.2	48	4.41
Bupropion	240.1	184	4	4.80	Normeperidine	234.1	56.1	20	4.41
Bupropion	240.1	131	20	4.80	Norquetiapine	296.1	210.1	24	5.38
Butocarboxim	213	156	5	5.14	Norquetiapine	296.1	139.1	60	5.38

化合物名称	母离子 (m/z)	子离子 (m/z)	碰撞能量 (V)	保留时间 (min)	化合物名称	母离子 (m/z)	子离子 (m/z)	碰撞能量 (V)	保留时间 (min)
Butocarboxim	213	75	10	5.14	Norsertaline	275	159.1	16	6.33
Caffeine	195.1	138	20	3.20	Norsertaline	275	89	72	6.33
Caffeine	195.1	83	40	3.20	Norverapamil	441.2	165	20	5.94
Cambendazole	303.1	261	16	5.57	Norverapamil	441.2	150.3	36	5.94
Cambendazole	303.1	217.1	32	5.57	Olaquinox	264	212	20	2.43
Carazolol	299.2	222	15	4.74	Olaquinox	264	143	30	2.43
Carazolol	299.2	116	20	4.74	Omeprazole	346.1	198.1	4	5.46
Carbadox	263	231	5	3.54	Omeprazole	346.1	136	28	5.46
Carbadox	263	90	30	3.54	Orbifloxacin	396.1	352.1	15	3.95
Carbamazepine	237	194	15	5.84	Orbifloxacin	396.1	295.1	22	3.95
Carbamazepine	237	179	35	5.84	Oxacillin	402.1	243.1	8	5.12
Carbamazepine 10,11 epoxide	253.1	210.2	8	4.96	Oxacillin	402.1	160	12	5.12
Carbamazepine 10,11 epoxide	253.1	180.1	24	4.96	Oxamyl	237	120	10	5.36
Carbaryl	202	145	5	6.38	Oxamyl	237	118	10	5.36
Carbaryl	202	117	10	6.38	Oxazepam	287	268.9	8	5.97
Carbofuran	222	165.1	10	6.21	Oxazepam	287	240.9	16	5.97
Carbofuran	222	123.1	20	6.21	Oxcarbazepine	253.1	208	16	5.25
Carbosulfan	381.1	160	15	11.53	Oxcarbazepine	253.1	180.1	24	5.25
Carbosulfan	381.1	118	15	11.53	Oxfendazole	316.1	191	17	4.92
Carisoprodol	261.2	176.2	0	6.20	Oxfendazole	316.1	159	35	4.92
Carisoprodol	261.2	55.1	24	6.20	Oxfendazole-Sulfone/ Fenbendazole sulfone	332.1	300	19	5.61
Carnidazole	245.1	118	10	5.01	Oxfendazole-Sulfone/ Fenbendazole sulfone	332.1	159	40	5.61
Carnidazole	245.1	75	40	5.01	Oxibendazole	250.1	218.1	14	5.92
Cetirizine	389.2	201	20	5.95	Oxibendazole	250.1	176	27	5.92
Cetirizine	389.2	166.1	40	5.95	Oxolinic acid	262.1	216	30	5.02
Chlormadinone acetate	405.2	345.2	8	8.83	Oxolinic acid	262.1	160	40	5.02
Chlormadinone acetate	405.2	309.2	12	8.83	Oxycodone	316.1	298.2	8	3.28
Chlorpheniramine	275.1	230.1	20	5.06	Oxycodone	316.1	241.1	24	3.28
Chlorpheniramine	275.1	167.1	40	5.06	Oxymorphone	302.1	284.1	12	2.27
Chlorpromazine	319.2	246	20	6.51	Oxymorphone	302.1	227	20	2.27
Chlorpromazine	319.2	86	15	6.51	Oxymorphone glucuronide	478.1	284.1	28	0.75
Chlorpyrifos	350	198	15	10.35	Oxymorphone glucuronide	478.1	227.1	48	0.75
Chlorpyrifos	350	79	35	10.35	Paracetamol/ 4-Acetamidophenol	152.1	110	13	2.52
Cinoxacin	263	217	20	4.24	Paracetamol/ 4-Acetamidophenol	152.1	65	32	2.52
Cinoxacin	263	189	30	4.24	Paraxanthine	181.1	124.1	20	2.67
Ciprofloxacin	332.1	314.1	20	3.67	Paraxanthine	181.1	42	40	2.67
Clenbuterol	277.1	203	12	4.16	Paroxetine	330.1	192.1	16	5.71
Clindamycin	425.2	377	20	4.95	Paroxetine	330.1	70.1	32	5.71
Clindamycin	425.2	126	30	4.95	Penicilline G	335.1	176	6	2.91
Clobetasol 17-propionate	467.2	372.9	6	8.55	Penicilline G	335.1	160	5	2.91

化合物名称	母离子 (m/z)	子离子 (m/z)	碰撞能量 (V)	保留时间 (min)	化合物名称	母离子 (m/z)	子离子 (m/z)	碰撞能量 (V)	保留时间 (min)
Clobetasol 17-propionate	467.2	354.9	8	8.55	Phenacetin	180	138	11	4.99
Clobetasone 17-butyrate	479.2	342.8	12	9.26	Phenacetin	180	110	19	4.99
Clobetasone 17-butyrate	479.2	278.9	14	9.26	Phenazone/Antipyrine	189	147	20	3.98
Clopidogrel carboxylic acid	308	198.1	8	4.21	Phenazone/Antipyrine	189	77	30	3.98
Clopidogrel carboxylic acid	308	76.9	64	4.21	Phenmetrazine	178.1	115	32	3.33
Clopidol	192	101	30	2.95	Phenmetrazine	178.1	91	36	3.33
Clopidol	192	87	35	2.95	Phentermine	150.1	133.1	4	3.51
Cloxacillin	436	277	10	5.11	Phenylbutazone	309	160	20	7.76
Cloxacillin	436	160	6	5.11	Phenylbutazone	309	77	40	7.76
Cocaethylene	318.2	196.2	12	4.90	Phenylpropanolamine	152.11	134	4	2.55
Cocaethylene	318.2	82.2	28	4.90	Phenylpropanolamine	152.11	117	12	2.55
Cocaine	304.1	182.2	16	4.47	Pioglitazone	357.13	133.9	24	7.16
Cocaine	304.1	77	64	4.47	Pioglitazone	357.13	119	48	7.16
Codeine	300.2	165.1	40	2.99	Pipracil	518.2	160	15	3.94
Codeine	300.2	128.1	60	2.99	Pipracil	518.2	143	25	3.94
Cortisone	361.2	163.1	20	5.46	Pirimicarb	239	182	15	6.41
Cortisone	361.2	121	30	5.46	Pirimicarb	239	72	20	6.41
Cotinine	177.1	98.1	24	2.99	Prednicarbate	489.2	380.9	6	8.43
Cotinine	177.1	80.1	28	2.99	Prednicarbate	489.2	114.8	12	8.43
CV	372.2	356.2	42	8.17	Prednisolone	361.2	343.1	6	5.33
CV	372.2	340.2	55	8.17	Prednisolone	361.2	146.9	20	5.33
Cyproheptadine	288.3	191	28	6.10	Prednisone	359.2	341.1	6	5.35
Cyproheptadine	288.3	96	24	6.10	Prednisone	359.2	147	24	5.35
Dapsone	249.1	156	9	4.37	Pregabalin	160.1	142.2	8	2.35
Dapsone	249.1	92	23	4.37	Pregabalin	160.1	55.2	20	2.35
DEET	192.1	119.3	16	6.58	Primidone	219.1	162.1	4	3.95
DEET	192.1	91	32	6.58	Primidone	219.1	91.2	24	3.95
Deflazacort	442.2	141.9	36	6.73	Progesterone	315.2	109	25	8.75
Deflazacort	442.2	123.9	50	6.73	Progesterone	315.2	97	21	8.75
Dehydroaripiprazole	446.1	285.1	24	6.33	Promethazine	285.1	86.1	16	5.82
Dehydroaripiprazole	446.1	98.1	44	6.33	Promethazine	285.1	71.1	48	5.82
Demeton	281	89	24	7.07	Propamocarb	189	144	10	2.73
Demeton	259	89	14	7.07	Propamocarb	189	102	15	2.73
Desloratadine	311.1	259.1	20	5.35	Propionylpromazine	341.2	86	20	6.25
Desloratadine	311.1	258.1	40	5.35	Propionylpromazine	341.2	58	30	6.25
Desmethylcitalopram	311.1	262.2	8	5.37	Propoxur	210	168	5	6.13
Desmethylcitalopram	311.1	109.1	20	5.37	Propoxur	210	111	10	6.13
Desmethylvenlafaxine	264.2	246.2	4	4.13	Propranolol	260.2	116.1	12	5.00
Desmethylvenlafaxine	264.2	58.1	16	4.13	Propranolol	260.2	56.1	24	5.00
Dexamethasone	393.1	355	4	5.95	Propyl thiouracil	171	154	14	3.40
Dexamethasone	393.1	146.8	25	5.95	Propyl thiouracil	171	112	17	3.40
Dextromethorphan	272.2	171.1	36	5.19	Pseudoephedrine	166.1	115.1	24	2.92
Dextromethorphan	272.2	128.1	64	5.19	Pseudoephedrine	166.1	91	32	2.92

化合物名称	母离子 (m/z)	子离子 (m/z)	碰撞能量 (V)	保留时间 (min)	化合物名称	母离子 (m/z)	子离子 (m/z)	碰撞能量 (V)	保留时间 (min)
Dichlorvos	221	145	15	4.07	Quetiapine	384.2	253.1	16	5.75
Dichlorvos	221	109	15	4.07	Quetiapine	384.2	221	36	5.75
Dicloxacillin	470	310.9	8	5.34	Quinoxaline-2-carboxylic Acid	175	129	13	2.32
Dicloxacillin	470	160	8	5.34	Quinoxaline-2-carboxylic Acid	175	102	33	2.32
Diethofencarb	268	226	5	7.60	Ractopamine	302	164.1	10	3.70
Diethofencarb	268	152	20	7.60	Ractopamine	302	121	22	3.70
Diflorasone diacetate	495.2	316.8	8	7.79	Rafoxanide	625.8	373	32	10.15
Diflorasone diacetate	495.2	278.8	10	7.79	Rafoxanide	625.8	127	40	10.15
Diltiazem	415.1	178.1	20	5.64	Ritalinic acid	220.1	84.2	16	3.36
Diltiazem	415.1	109.1	76	5.64	Ritalinic acid	220.1	56.1	40	3.36
Dimethoate	230	199	5	4.60	Ronidazole	201.1	140	4	2.94
Dimethoate	230	171	10	4.60	Ronidazole	201.1	55	18	2.94
Dimetridazole	142.1	96	14	3.41	Salbutamol	240	222.1	5	2.39
Dimetridazole	142.1	81	30	3.41	Salbutamol	240	148	13	2.39
Dioxacarb	224	167	5	4.66	Salinomycin	768.5	733.4	15	12.43
Dioxacarb	224	123	15	4.66	Salinomycin	768.5	715.5	20	12.43
Diphenhydramine	256.2	167.1	4	5.44	Secnidazole	186.1	128.1	15	3.51
Diphenhydramine	256.2	165.1	44	5.44	Secnidazole	186.1	59	20	3.51
Dipyron	310.1	191	4	7.78	Sertraline	306.1	275	4	6.34
Dipyron	310.1	80	32	7.78	Sertraline	306.1	159.1	20	6.34
Disopyramide	340.2	239.1	8	4.38	Sildenafil	475.2	100	24	6.09
Disopyramide	340.2	194.2	48	4.38	Sildenafil	475.2	58.1	40	6.09
Donepezil	380.2	91.1	40	5.11	Simvastatin	419.3	198.9	12	9.84
Donepezil	380.2	65.2	76	5.11	Simvastatin	419.3	173	28	9.84
Doramectin	921.4	777.4	45	11.62	Sineptina-772	772.5	174	35	6.45
Doramectin	921.4	449.2	55	11.62	Sineptina-772	772.5	109	45	6.45
Doxepin	280.2	107.1	20	5.58	Sineptina-786	786.5	229.3	45	6.96
Doxepin	280.2	77.1	60	5.58	Sineptina-786	786.5	109	45	6.96
Duloxetine	298.1	153.9	0	5.93	Sotalol	273.1	255.1	4	2.55
Duloxetine	298.1	44	12	5.93	Sotalol	273.1	133.1	20	2.55
Ecgonine methyl ester	200.1	182	12	0.77	Spiramycin	843.5	174	42	5.43
Ecgonine methyl ester	200.1	82.1	24	0.77	Spiramycin	843.5	101	46	5.43
Econazole	381	125	40	9.25	Sulfabenzamide	277.1	156	10	4.35
Econazole	381	69	20	9.25	Sulfabenzamide	277.1	108	25	4.35
EDDP	278.2	249	20	5.78	Sulfachloropyridazine	285	156	10	4.38
EDDP	278.2	234.1	24	5.78	Sulfachloropyridazine	285	108	25	4.38
Epitestosterone	289.2	109	22	7.56	Sulfadiazine	251.1	156	10	3.07
Epitestosterone	289.2	97	24	7.56	Sulfadiazine	251.1	108	22	3.07
Eprinomectin	936.4	490.3	59	10.45	Sulfadimethoxine	311	156	20	5.31
Eprinomectin	936.4	352.2	65	10.45	Sulfadimethoxine	311	108	26	5.31
Erythromycin	734.5	576.3	14	5.46	Sulfadimidine	279.1	186.1	15	4.01
Erythromycin	734.5	158.1	30	5.46	Sulfadimidine	279.1	156.1	16	4.01
Erythromycin-anhydro	716.4	158	24	5.77	Sulfadoxine	311.1	156	15	4.65

化合物名称	母离子 (m/z)	子离子 (m/z)	碰撞能量 (V)	保留时间 (min)	化合物名称	母离子 (m/z)	子离子 (m/z)	碰撞能量 (V)	保留时间 (min)
Erythromycin-anhydro	716.4	83.2	76	5.77	Sulfadoxine	311.1	92	30	4.65
Escitalopram	325.2	262.2	16	5.46	Sulfaguanidine	215	156	10	1.33
Escitalopram	325.2	109.1	20	5.46	Sulfaguanidine	215	108	20	1.33
Famotidine	338	189.1	12	2.51	Sulfamerazine	265.1	172	12	3.61
Famotidine	338	155	28	2.51	Sulfamerazine	265.1	156	15	3.61
Fenbendazole	300.1	268	18	7.22	Sulfamerter/ Sulfamethoxy pyridazine	281	156	15	3.99
Fenbendazole	300.1	159	36	7.22	Sulfamerter/ Sulfamethoxy pyridazine	281	108	25	3.99
Fenobucarb	208.2	152.1	5	7.44	Sulfamethazine	279.1	186	8	3.97
Fenobucarb	208.2	95	10	7.44	Sulfamethazine	279.1	92.1	28	3.97
Fenoterol	304.1	286.2	8	2.97	Sulfamethizole	271	156	10	3.73
Fenoterol	304.1	135.2	15	2.97	Sulfamethizole	271	108	22	3.73
Fenoxycarb	302.1	116	10	8.27	Sulfamethoxazole	254.1	156	10	4.64
Fenoxycarb	302.1	88	20	8.27	Sulfamethoxazole	254.1	108	25	4.64
Fentanyl	337.2	188.3	20	5.45	Sulfamonomethoxine	281.1	156.1	15	4.27
Fentanyl	337.2	105.1	36	5.45	Sulfamonomethoxine	281.1	108.1	26	4.27
Flubendazole	314.1	282.1	18	6.23	Sulfamoxol	268	156	13	4.52
Flubendazole	314.1	123	37	6.23	Sulfamoxol	268	113	16	4.52
Fluconazole(l) (Diflucan)	307.1	238	10	4.11	Sulfanitran	336.1	293.9	8	6.07
Fluconazole(l) (Diflucan)	307.1	220	20	4.11	Sulfanitran	336.1	156	8	6.07
Fludrocortisone 21-acetate	423.2	238.9	22	6.62	Sulfaphenazole	315	222	15	5.38
Fludrocortisone 21-acetate	423.2	120.9	36	6.62	Sulfaphenazole	315	158	30	5.38
Fludroxycortide/Flurandrenolide	437.2	180.9	30	6.47	Sulfapyrazole	329.1	172	30	5.65
Fludroxycortide/Flurandrenolide	437.2	120.8	40	6.47	Sulfapyrazole	329.1	145	50	5.65
Flumequine	262	202	30	6.07	Sulfapyridine	250.1	184	15	3.45
Flumequine	262	126	50	6.07	Sulfapyridine	250.1	156	10	3.45
Flumethasone	411.2	253	10	6.00	Sulfaquinoxaline	301.1	156	11	5.26
Flumethasone	411.2	121.1	34	6.00	Sulfaquinoxaline	301.1	108	22	5.26
Flumethasone pivalate	495.3	475.2	10	8.66	Sulfathiazole	256	156	10	3.28
Flumethasone pivalate	495.3	57.1	20	8.66	Sulfathiazole	256	108	21	3.28
Flunixin/Flunixin Meglumine	297.1	279.1	32	6.27	Sulfisomidine	279	186	15	2.88
Flunixin/Flunixin Meglumine	297.1	264.1	4	6.27	Sulfisomidine	279	124	20	2.88
Fluocinolone acetonide	453.2	413.2	10	6.40	Sulphacetamide	215.1	156.1	5	2.60
Fluocinolone acetonide	453.2	121.1	40	6.40	Sulphacetamide	215.1	92	20	2.60
Fluoromethalone	377.2	320.9	8	6.59	Sulphisoxazole/Sulfisoxazole/ Sulfafurazole	268.1	156	10	3.78
Fluoromethalone	377.2	278.9	10	6.59	Sulphisoxazole/Sulfisoxazole/ Sulfafurazole	268.1	113	10	3.78
Fluoxetine	310.1	148.1	2	6.28	Sulpiride	342.1	214	40	2.74
Fluoxetine	310.1	44	8	6.28	Sulpiride	342.1	112.1	20	2.74
Fluphenazine	438.1	171.1	26	6.95	Sumatriptan	296.1	155.9	52	3.12
Fluphenazine	438.1	143	30	6.95	Sumatriptan	296.1	58	12	3.12
Fluticasone propionate	501.2	312.9	8	8.60	Tadalafil	390.1	268	0	6.32

化合物名称	母离子 (m/z)	子离子 (m/z)	碰撞能量 (V)	保留时间 (min)	化合物名称	母离子 (m/z)	子离子 (m/z)	碰撞能量 (V)	保留时间 (min)
Fluticasone propionate	501.2	292.9	10	8.60	Tadalafil	390.1	204	80	6.32
Formoterol	345	149	15	4.10	Temazepam	301.1	283	4	6.67
Formoterol	345	121	20	4.10	Temazepam	301.1	255	16	6.67
Furathiocarb	383	252	10	9.92	Terbutaline	226.1	152.1	10	2.35
Furathiocarb	383	195	15	9.92	Terbutaline	226.1	107	25	2.35
Gabapentin	172.1	154.1	8	2.37	Terfenadine	472.3	454.3	20	7.55
Gabapentin	172.1	55	24	2.37	Terfenadine	472.3	436.3	20	7.55
Glyburide	494.1	369	12	7.71	Testosterone	289.2	109	23	7.02
Glyburide	494.1	169.1	36	7.71	Testosterone	289.2	97	21	7.02
Griseofulvin	353.1	215	20	6.73	Thiabendazole	202	175	25	4.72
Griseofulvin	353.1	165.1	20	6.73	Thiabendazole	202	131	35	4.72
Halcinonide	455.2	121	40	8.42	Thiodicarb	355.1	163.1	12	6.01
Halcinonide	455.2	104.9	48	8.42	Thiodicarb	355.1	88.1	27	6.01
Halofuginone	414	138.1	16	4.84	Thiofanox	241	184	5	6.49
Halofuginone	414	100.1	16	4.84	Thiofanox	241	57.1	20	6.49
Haloperidol	376.1	165.1	20	5.74	Tilmicosin	869.6	696.4	45	6.15
Haloperidol	376.1	123	40	5.74	Tilmicosin	869.6	174	50	6.15
HMMNI/2-Hydroxymethyl-1-methyl-5-nitroimidazole	158.1	140	10	2.64	Tinidazole	248	121	15	3.70
HMMNI/2-Hydroxymethyl-1-methyl-5-nitroimidazole	158.1	55	18	2.64	Tinidazole	248	93	20	3.70
Hydrocodone	300.1	199.1	28	3.41	Tramadol	264.2	58.1	12	4.13
Hydrocodone	300.1	171.1	40	3.41	Tramadol	264.2	56.1	68	4.13
Hydrocortisone	363.2	121	24	5.40	Trazadone	372.1	176.1	20	5.45
Hydrocortisone	363.2	105.1	50	5.40	Trazadone	372.1	148.1	36	5.45
Hydromorphone	286.1	185.1	24	2.52	Triamcinolone	395.2	357.1	8	4.22
Hydromorphone	286.1	157.1	48	2.52	Triamcinolone	395.2	225.1	14	4.22
Hydroxybupropion	256.1	238.1	4	4.15	Triamcinolone acetonide	435.2	396.9	10	6.27
Hydroxybupropion	256.1	130.1	48	4.15	Triamcinolone acetonide	435.2	338.9	10	6.27
Hydroxyzine	375.2	201	20	6.14	Triamterene	254.1	237.1	24	3.66
Hydroxyzine	375.2	166.1	40	6.14	Triamterene	254.1	104.1	32	3.66
Imipramine	281.2	86.1	12	6.05	Triclabendazole	359	343.9	24	8.79
Imipramine	281.2	58.2	44	6.05	Triclabendazole	359	274	40	8.79
Indoxacarb	528	293	10	9.37	Trimethoprim	291.1	230.1	25	3.50
Indoxacarb	528	249	10	9.37	Trimethoprim	291.1	123	25	3.50
Iprnidazole	170.1	124	16	5.12	Tripelenamine	256.2	211.1	10	5.21
Iprnidazole	170.1	109	26	5.12	Tripelenamine	256.2	91.1	40	5.21
Iprnidazole-OH/Hydroxy-Iprnidazole	186.1	168	10	4.02	Tulobuterol	228.1	172	5	4.12
Iprnidazole-OH/Hydroxy-Iprnidazole	186.1	122	20	4.02	Tulobuterol	228.1	154	13	4.12
Isoprocarb	194	137	5	6.83	Tylosin	916.5	174	42	5.83
Isoprocarb	194	95	20	6.83	Tylosin	916.5	101	54	5.83
Ivermectin	897.4	753.3	46	12.26	Valsartan	436.2	235.1	12	5.85

化合物名称	母离子 (m/z)	子离子 (m/z)	碰撞能量 (V)	保留时间 (min)	化合物名称	母离子 (m/z)	子离子 (m/z)	碰撞能量 (V)	保留时间 (min)
Ivermectin	897.4	329.2	60	12.26	Valsartan	436.2	207.2	20	5.85
Ketoconazole	531.2	489.1	40	7.36	Venlafaxine	278.2	260.3	0	4.69
Ketoconazole	531.2	82.1	40	7.36	Venlafaxine	278.2	58.2	20	4.69
Ketoprofen	255.1	209.2	4	5.88	Verapamil	455.3	165.1	24	6.07
Ketoprofen	255.1	77.1	52	5.88	Verapamil	455.3	150.1	40	6.07
Lamotrigine	256	109	52	4.25	Virginiamycin M1	526.3	508.4	15	6.53
Lamotrigine	256	74	76	4.25	Virginiamycin M1	526.3	355	20	6.53
LCV	374.3	358.2	30	11.01	Xylazine	221.1	164	24	4.14
LCV	374.3	238.1	26	11.01	Xylazine	221.1	90	21	4.14
Levamisole	205	178	20	3.20	Zolpidem	308.2	236.1	24	5.51
Levamisole	205	91	30	3.20	Zolpidem	308.2	235.1	36	5.51
Levorphanol	258.2	199.1	20	3.95	Zolpidem phenyl-4-carboxylic acid	338.1	265	36	3.48
Levorphanol	258.2	157.2	32	3.95	Zolpidem phenyl-4-carboxylic acid	338.1	65.1	76	3.48

表 5. 正离子模式下化合物 MRM 离子对参数列表

化合物名称	母离子 (m/z)	子离子 (m/z)	碰撞能量 (V)	保留时间 (min)	化合物名称	母离子 (m/z)	子离子 (m/z)	碰撞能量 (V)	保留时间 (min)
11-nor-9-carboxy-delta-THC	343.2	299.3	20	6.10	Furosemide	329	285	8	4.26
11-nor-9-carboxy-delta-THC	343.2	245	24	6.10	Furosemide	329	204.9	16	4.26
17alpha-Estradiol	271.1	183	45	4.75	Gemfibrozil	249.1	127.1	4	5.85
17alpha-Estradiol	271.1	145	45	4.75	Gemfibrozil	249.1	121	4	5.85
17beta-Estradiol	271.1	183	45	4.9	Hexestrol	269.1	133	15	5.18
17beta-Estradiol	271.1	145	45	4.9	Hexestrol	269.1	119	43	5.18
2,4,5-T	255	197	10	4.91	Hydrochlorothiazide	295.9	268.9	12	2.89
2,4,5-T	253	195	10	4.91	Hydrochlorothiazide	295.9	205.2	20	2.89
2,4,5-TP	269	197	8	5.23	Ibuprofen	205	161	2	5.49
2,4,5-TP	267	195	8	5.23	MCPA	201	143	10	4.51
2,4-D	221	163	15	4.45	MCPA	199	141	10	4.51
2,4-D	219	161	15	4.45	MCPB	229	143	10	5.20
2,4-DB	249	163	10	5.18	MCPB	227	141	10	5.20
2,4-DB	247	161	10	5.18	MCPB	215	143	10	4.88
2,4-DP	235	163	5	4.82	MCPB	213	141	10	4.88
2,4-DP	233	161	5	4.82	MCPB	151	135.9	8	3.76
Bezafibrate	360.1	274	12	4.81	Methylparaben	151	92	16	3.76
Bezafibrate	360.1	154.1	28	4.81	Modafinil acid	273	167.1	8	4.17
Celecoxib	380.1	316.1	20	5.50	Modafinil acid	273	165	36	4.17
Celecoxib	380.1	276	28	5.50	Naproxen	229.1	170.1	4	4.78
Chloramphenicol	321	257	10	3.77	Naproxen	229.1	168.9	28	4.78
Chloramphenicol	321	152	15	3.77	n-Butylparaben	193.1	136	12	5.01
Diclozauril	407	336	15	5.53	n-Butylparaben	193.1	92	20	5.01
Diclozauril	405	334	14	5.53	Nicarbazin	301	137	5	5.23
Diclofenac	294	250	8	5.41	Nicarbazin	301	107	35	5.23
Diclofenac	294	35	32	5.41	Nitroxynil	288.9	162	20	5.02
Diclofenac 4-hydroxy	310	265.9	8	4.62	Nitroxynil	288.9	127	28	5.02
Diclofenac 4-hydroxy	310	34.7	32	4.62	Phenobarbital	231.1	188.1	0	3.73
Dienestrol	265.1	249	25	5.17	Phenobarbital	231.1	42.1	12	3.73
Dienestrol	265.1	93	30	5.17	Phenytoin	251.1	102.1	20	4.15
Diethylstilbestrol	267.1	251	27	5.11	Phenytoin	251.1	41.7	60	4.15
Diethylstilbestrol	267.1	237	30	5.11	Pravastatin	423.2	320.9	12	3.88
Estriol	287.1	171	43	3.66	Pravastatin	423.2	303.2	16	3.88
Estriol	287.1	145	50	3.66	Thiamphenicol	354	290	6	3.11
Ethynyl estradiol	295.2	269	35	4.95	Thiamphenicol	354	185	20	3.11
Ethynyl estradiol	295.2	145	45	4.95	Triclocarban	312.9	160.1	8	6.05
Fenbufen	253.1	153.1	20	4.87	Triclocarban	312.9	35	44	6.05
Fenbufen	253.1	55	24	4.87	Triclosan	286.9	35	4	6.07
Florfenicol	356	336	5	3.68	Warfarin	307.1	250	16	5.08
Florfenicol	356	185	15	3.68	Warfarin	307.1	161.2	16	5.08

结果与讨论

通过分析混合标准溶液对仪器方法进行优化,可在 20 min 内对 377 种 PPCP 进行有效分离。各化合物响应良好,绝大多数化合物检测限可低于 10 ng/L,灵敏度足够高,无需进行样品富集。图 1 显示了 100 ng/L PPCP 混合标准样品的提取离子色谱叠加图,保留时间如表 4 和表 5 所示。

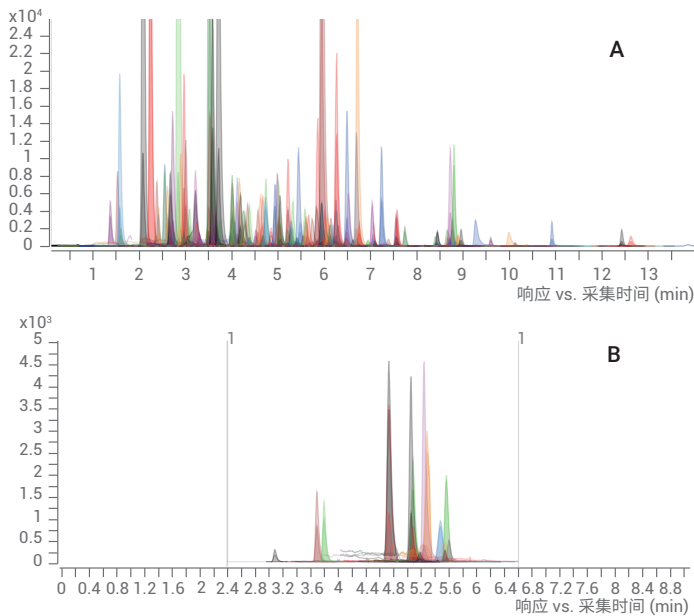


图 1. 377 种 PPCP 混合标准样溶液 (浓度 100 ng/L) 在正离子模式 (A) 和负离子模式 (B) 下的提取离子色谱叠加图

采用本文所述方法在浓度范围 10 ng/L 到 10^5 ng/L 内对 377 种 PPCP 的线性相关性进行讨论。如图 2 所示,在正离子模式下阿特拉津和负离子模式下布洛芬的校准曲线的线性相关系数 (R^2) 分别为 0.996 和 0.995,线性关系良好,大多数化合物的线性相关系数均大于 0.99。

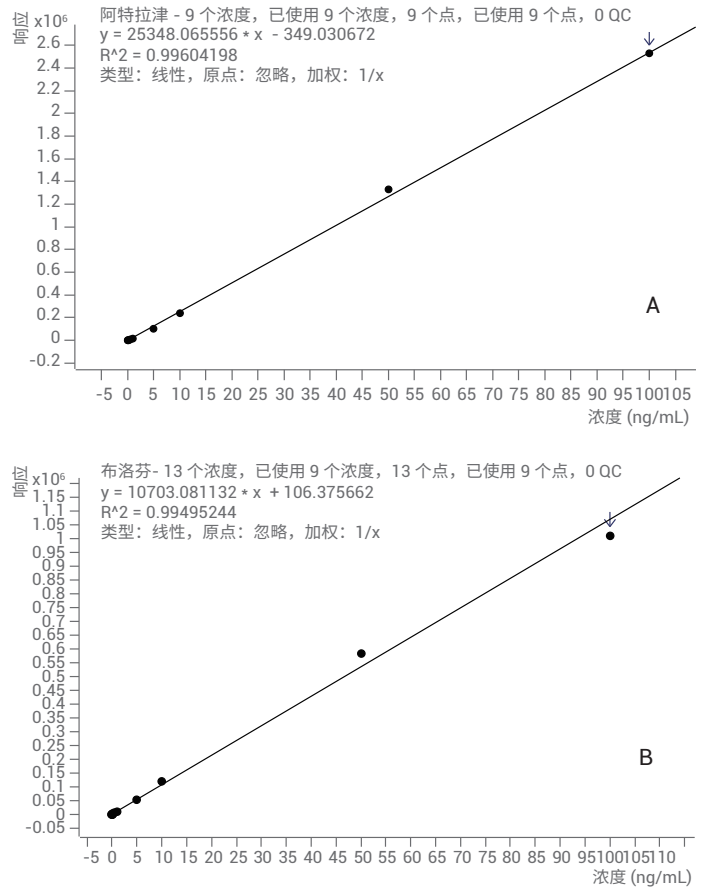


图 2. 阿特拉津 (A) 和布洛芬 (B) 的校准曲线 (浓度范围 10 - 10^5 ng/L)

浓度为 100 ng/L 实际水样加标样品,连续进样 7 针,计算各化合物的相对标准偏差。如图 3 所示,超过 78% 的化合物相对 RSD 小于 20%,具有良好的重复性。

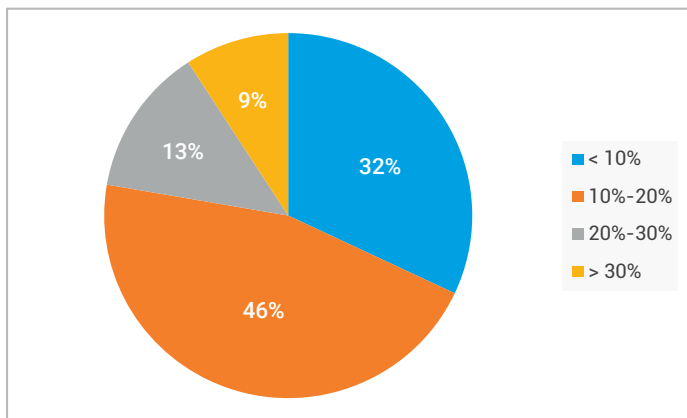


图 3. 100 ng/L 添加浓度下连续进样 7 针所得各化合物 RSD 分布图

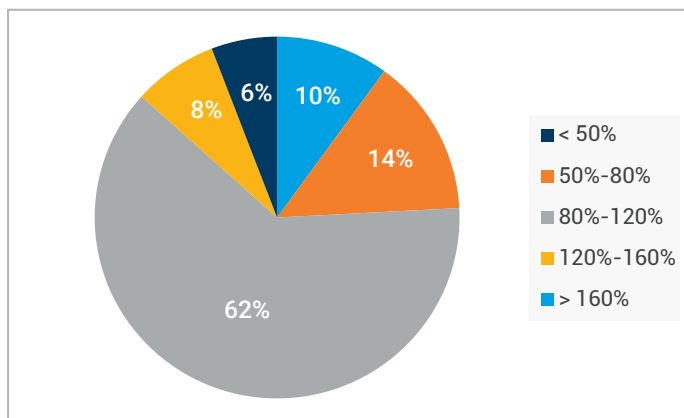


图 4. 100 ng/L 添加浓度下各化合物回收率分布图

浓度为 100 ng/L 实际水样加标样品，按方法进样定量分析，计算回收率。如图 4 所示，大多数样品回收率在 80%-120% 之间，回收率良好。

实际水样取自北京某湖泊及周边水域，在 10 个不同位置分别取 10 个水样，采用 0.22 μm 滤膜过滤后直接进样 20 μL 进行分析检测。如表 6 所示，共检测到 PPCP 共 13 个，其中尼泊金甲酯浓度超过 4 μg/L，林可霉素、舒必利、副黄嘌呤、阿特拉津、氟康唑、氟苯尼考部分水样中浓度超过 100 ng/L。

表 6. 实际水样 PPCP 检测结果（浓度单位：ng/L）

样品编号	扑热息痛	2-氨基苯并咪唑	副黄嘌呤	舒必利	可铁宁	林可霉素	氟康唑	磺胺甲恶唑	氯苯那敏	莠去津	益康唑	尼泊金甲酯	氟甲砜霉素
1	-	5.8	45.8	47.0	19.4	82.6	43.0	24.5	13.5	84.0	43.4	9336.2	68.2
2	-	4.7	45.3	11.5	10.7	31.9	64.3	17.8	12.9	70.6	14.2	4548.3	31.0
3	4.3	-	-	3.6	8.3	10	56.5	10.0	18.1	46.9	7.3	8265.4	9.1
4	4.9	2.7	38.4	23.8	19.5	51.9	144	17.1	11.9	24.0	8.7	9906.1	15.6
5	-	9.5	29.5	53.8	28.8	104	69.6	49.1	10.6	54.6	14.7	4289.8	127.4
6	1.7	3.8	43.8	22.8	20.5	66.8	53.1	-	11.5	33.2	22.6	9879.2	12.6
7	-	7.1	69.1	90.2	23.8	160	79.9	40.7	14.2	62.6	5.4	7010.9	279.5
8	92.7	8.3	144.7	88.4	25.7	164.3	85.9	39.8	40.1	55.9	10.0	6581.0	146.9
9	-	6.1	42.5	52.7	20.3	99.8	64.2	33.9	11.2	31.3	8.9	5288.0	35.5
10	4.5	21.2	285.5	137.4	6.6	560.7	70.7	54.1	40.1	240.8	5.7	5420.9	197.9

结论

本文采用 Agilent 1290 Infinity II 液相色谱/6495 三重四极杆液质联用系统，应用动态 MRM 方式成功实现了同时对环境水中 377 种 PPCP 的高灵敏度、快速、简单、准确的筛查和定量分析。本方法通过直接进样方法即可满足常规检测要求，有助于省掉繁琐的样品富集和净化步骤，从而显著提高样品分析通量。

查找当地的安捷伦客户中心：

www.agilent.com/chem/contactus-cn

免费专线：

800-820-3278, 400-820-3278 (手机用户)

联系我们：

LSCA-China_800@agilent.com

在线询价：

www.agilent.com/chem/erfq-cn

www.agilent.com

安捷伦对本资料可能存在的错误或由于提供、展示或使用本资料所造成的间接损失不承担任何责任。

本文中的信息、说明和技术指标如有变更，恕不另行通知。

© 安捷伦科技（中国）有限公司，2017
2017年10月22日，中国印刷
5991-8660ZHCN

