



# Analysis of Pesticide Residues in Mango by GC/MS/MS With Bond Elut QuEChERS EN Kits

## Application Note

Food Testing and Agriculture

### Authors

Claudia Piscato, Claudia Barbosa, and  
Amir Gebara

Instituto Biológico  
Laboratório de Resíduos de Pesticidas  
São Paulo, SP, Brazil

Kumi Shiota Ozawa  
Agilent Technologies Brasil Ltda.  
Barueri, SP, Brazil

### Abstract

This application note describes the quantitative analysis of 19 of the 40 regulated pesticides with maximum residue limits (MRLs) established by the Brazilian Health Surveillance Agency (ANVISA). We also analyzed nine pesticides found in mango during the monitoring programs developed in the Pesticide Residues Laboratory (LRP/IB)/São Paulo State and by governmental programs. The samples were also analyzed with a qualitative multiresidue method for 258 pesticides, extending the analytical scope to analytes with low probability of being present, as recommended in SANCO/12571/2013 guidelines [1]. The extraction was performed using an Agilent Bond Elut QuEChERS EN kit. Target pesticides were analyzed by GC/MS/MS using an Agilent 7890A GC and an Agilent 7000B Triple Quadrupole GC/MS in a constant pressure/postcolumn backflush configuration (Pesticide Analyzer 411). The method was validated in terms of recovery and reproducibility. The limits of detection (LOD) ranged between 0.0006 and 0.0607 mg/kg and limits of quantitation (LOQ) were between 0.0025 and 0.5 mg/kg. LOQs were established as  $\frac{1}{4}$  of MRL. Recoveries for all compounds were 70 to 120%, and RSDs were below 20% for five replicates. LODs were calculated as three times the RSDs at LOQ levels.



**Agilent Technologies**

## Introduction

Mango is one of the most consumed fresh fruits in the world, and Brazil is one of the main exporters. It is a fruit with a large range of varieties (Palmer, Tommy Atkins, Edward, Peter, Haden, Keith, Mummy, Julie, Saigon, Dabsha, Hindi, Van Dyke, Benisha, Kesar, and Lick). Each one has its own characteristics in terms of pulp quality, taste, color, and so on. In general, mango is consumed fresh and unprocessed. However, it is possible to find it in different forms, such as purée, juice, sliced, chutney, and flakes. World market demand for mangoes has been rapidly growing. For example, growth averaged nine percent per year from 2006 to 2010 [2].

The international market varies considerably and depends on the preference of the consumer. To be exported, the fruit must exhibit a brilliant red color, have short fibers, and weigh around 250 to 600 g. In general, Tommy Atkins is the variety with a large market due mainly to its intense color and durability in long-distance transport [3]. For this reason, Tommy Atkins was chosen for this study.

Increasing the scope of the quantitative method was considered an illogical expense in terms of the cost of further standards and their application. However, to assess the presence of pesticides with low probability, the samples were also screened qualitatively with the same chromatographic method but with MRM transitions for 258 pesticides, including isomers and metabolites.

The Brazilian Health Surveillance Agency (ANVISA) regulates and lists maximum residue limits (MRLs) for about 40 pesticides for use on mango [4]. Of these regulated pesticides, 19 were selected on the basis of amenability to GC, and availability of standards. In addition, nine other pesticides found in mango during the national monitoring programs were included in the list, 28 compounds were analyzed quantitatively.

Two Brazilian national monitoring programs have begun, namely Pesticide Residues Analysis Program (PARA) on Food, developed by ANVISA, and National Plan for Control of Residues and Contaminants in Products of Plant Origin (PNCRC), developed by the Ministry of Agriculture. These initiatives are based on programs initially set up at the Instituto Biológico's Laboratório de Resíduos de Pesticidas (LRP) pesticide residues laboratory and São Paulo General Warehousing and Centers Company (CEAGESP) in 1978, with LRP participating in the PNCRC program. National programs were created to establish a service to evaluate and promote

food quality regarding the use of pesticides and similar compounds [4]. The high quantity of pesticides not allowed for specific crops was also verified in PARA and PNCRC monitoring programs, and from LRP analysis [4,9,10,11].

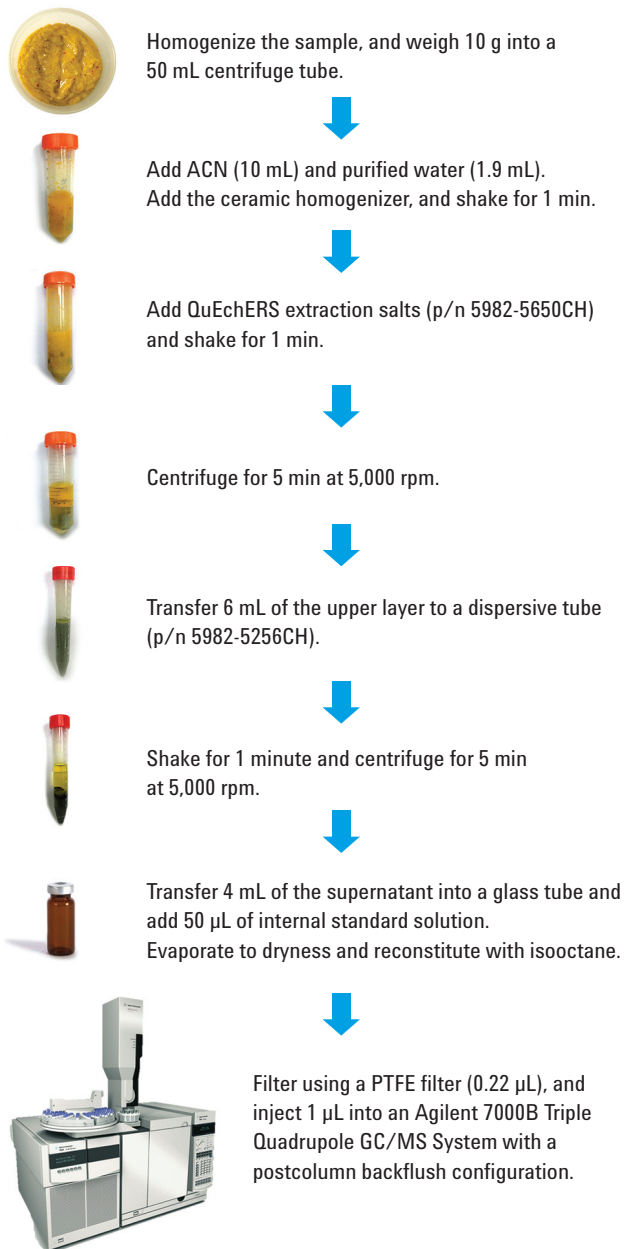
The programs list some commodities that may or may not be included in the annual sampling plan; mango is one of them. The others are rice, zucchini, pineapple, lettuce, banana, potato, beet, onion, carrot, cabbage, bean, orange, apple, papaya, corn, strawberry, cucumber, green pepper, cabbage, tomato, and grape [4].

## Materials and Methods

Acetonitrile, isooctane, and acetone were pesticide-residue grade. Pure standards from AccuStandard, around 99% pure, were used to prepare stock solutions at 1,000 ng/μL and working solutions that varied in concentration. Triphenyl phosphate at 0.5 mg/kg was prepared in isooctane and used as internal standard.

We performed extractions using the Agilent QuEChERS Extraction Kit for EN method 15662EN (p/n 5982-5650CH), in which 10 g of mango sample was extracted using premixed sachets of 4 g MgSO<sub>4</sub>, 1 g NaCl, 1 g Na citrate, and 0.5 g disodium citrate sesquihydrate. More details are shown in Figure 1. Mango is a highly pigmented fruit, so the GC components should be kept free from pigments and nonvolatiles [8]. We chose a subsequent dispersive cleanup designed to include pigment removal (Agilent Bond Elut QuEChERS SPE Dispersive Kit for Pigmented Fruits and Vegetables, EN method, p/n 5982-5256CH). This included premixed sachets containing 150 mg PSA, 15 mg graphitized carbon (GCB), and 885 mg MgSO<sub>4</sub>. Cleanup removed polar organic acids, some sugars and lipids, and carotenoids and chlorophyll. It is known that the use of GCB can affect planar pesticides such as quintozone and thiabendazole in our target list for quantitation [5]. However, the relatively low level of GCB in the Agilent formulation was a key criterion in its choice, and the good RSDs and recoveries we found (see Table 2) supported this approach.

A calibration in matrix extract was prepared daily and injected before and after the sample set to check calibration curve integrity. TPP was added after cleanup to avoid its retention by GCB [8]. The Agilent 7000B Triple Quadrupole GC/MS System was configured according to the Agilent Pesticide Analyzer 411 configuration, featuring a 30 m analytical column with postcolumn backflush.



## Instrumental conditions

### GC conditions

Column:	Agilent J&W DB-5ms, 30 m $\times$ 0.25 mm, 0.25 $\mu$ m (p/n 122-5532)
Inlet:	Split/splitless
Inlet liner:	Splitless, single taper, Ultra Inert liner with glass wool (p/n 5190-3167)
Carrier:	Helium
Inlet pressure:	36 psi (constant pressure mode) during run, 1 psi during backflush
Inlet temp:	280 $^{\circ}$ C
Inj vol:	1 $\mu$ L
Purge flow to split vent:	30 mL/min at 0.75 min
Gas saver:	On (20 mL/min at 2.0 min)
Oven temp :	70 $^{\circ}$ C (1 min), 50 $^{\circ}$ C/min to 150 $^{\circ}$ C (0 min), 6 $^{\circ}$ C/min to 200 $^{\circ}$ C (0 min), 16 $^{\circ}$ C/min to 280 $^{\circ}$ C (5.5 min)
Capillary flow technology:	Agilent Purged Ultimate Union (p/n G3186) used for backflushing the column and retention gap
Restrictor:	Deactivated capillary tubing, 0.7 m $\times$ 0.15 mm
Retention time locking:	Chlorpyrifos-methyl locked at 16.59 min
GC:	Agilent 7890A series (G3440A)
Autosampler:	Agilent 7693A Automatic Liquid Sampler injector and sample tray

### MS conditions

Spectrometer:	Agilent 7000B Triple Quadrupole GC/MS System
Mode:	Electron Impact
Transfer line temp:	280 $^{\circ}$ C
Solvent delay:	2.3 min
Source temp:	300 $^{\circ}$ C
Quadrupole temp:	Q1 and Q2 = 180 $^{\circ}$ C

Figure 1. Mango sample preparation process.

## Results and Discussion

### Performance evaluation, quantitative method

Blank samples were spiked with concentrations of 1 MRL, ½ MRL, and ¼ MRL. Calibration curves were prepared at five levels in concentrations of ¼, ½, 1, 2.5, and 5x MRL, all above  $R^2 = 0.99$  linearity. Recoveries ranged from 70 to 120% at ¼ MRL for the pesticides listed in Table 1.

In an effort to increase the low recovery of carbosulfan, 600 µL of 5 N NaOH were added after the addition of salts into the sample [6]. The addition of NaOH increased the pH from 4 to 6. Carbosulfan is an acid-sensitive pesticide; therefore, at pH 4 it may be hydrolyzed and converted into carbofuran. Carbosulfan exhibited increased recovery after pH adjustment from 59 to 113% when spiking at 0.0125 mg/kg. However, due to their sensitiveness to basic solutions [7], this addition affected the recoveries of vinclozolin, cyhalothrin, and cypermethrin, lowering them to about 50%. Therefore, it is recommended to prepare both solutions to have better recoveries for all pesticides.

Table 1. Pesticides analyzed with limits of detection (LOD), limits of quantitation (LOQ), RSD, recovery (REC), and MRM transitions.

No.	Pesticide	MRL	RT (min)	LOD	LOQ	RSD (%)	REC (%)	Quant	CE (V)	Qual	CE (V)
1	Azoxystrobin	0.3	36.6	0.0287	0.0750	13.1	97	344.1 → 329.0	15	344.1 → 182.9 344.1 → 171.9	25 40
2	BHC- <i>alpha</i>	0.01*	12.3	0.0007	0.0025	9.4	95	218.8 → 183.0	5	218.8 → 145.0 181.0 → 145.0	20 15
3	Bifenthrin	0.1	29.0	0.0088	0.0250	13.8	85	181.2 → 165.2	25	166.2 → 165.2 181.2 → 166.2	20 10
4	Carbosulfan	0.05	28.7	0.0016	0.0125	7.1	59/113*	118.0 → 76.0	5	164.0 → 103.1 164.0 → 149.0	25 10
5	Chlorpyrifos	0.05*	19.1	0.0047	0.0125	15.1	83	313.8 → 257.8	15	196.9 → 107.0	40
6	Cyhalothrin (λ)	0.1	30.5	0.0089	0.0250	12.6	94	208.0 → 181.0	7	197.0 → 161.0 197.0 → 141.0	5 10
7	Cypermethrin I	0.7	33.1	0.0480	0.1750	10.9	84	163.0 → 127.0	5	209.0 → 116.0 181.0 → 127.0 165.0 → 127.0	15 30 5
8	Diazinon	0.05*	14.4	0.0052	0.0125	12.4	112	137.1 → 84.0	10	137.1 → 54.0 199.1 → 93.0	20 15
9	Difenoconazole I	0.2	35.5	0.0170	0.0500	15.6	72	322.8 → 264.8	15	264.9 → 202.0 324.8 → 266.8	20 15
10	Endosulfan I	0.01*	23.1	0.0010	0.0025	13.1	100	239.0 → 204.0	15	339.0 → 267.0 207.0 → 172.0 241.0 → 136.0 241.0 → 206.0	2 15 20 20
11	Endosulfan II	0.01*	25.7	0.0010	0.0025	17.3	81	241.0 → 206.0	15	339.0 → 267.0 241.0 → 136.0 239.0 → 204.0 207.0 → 172.0	2 40 15 15
12	Endosulfan sulfate	0.01*	27.1	0.0011	0.0025	17.0	85	271.8 → 237.0	15	387.0 → 253.0 271.8 → 235.0	10 15
13	Ethion	0.05*	26.1	0.0049	0.0125	13.1	100	152.9 → 96.9	10	124.9 → 96.9 230.9 → 175.0	0 10
14	Etofenprox	0.3	33.5	0.0219	0.0750	12.7	77	163.0 → 107.1	20	135.0 → 107.0 163.0 → 135.1	10 10
15	Famoxadone	0.2	37.1	0.0163	0.0500	10.9	100	223.9 → 196.2	10	197.0 → 141.1 197.0 → 115.0	15 30
16	Fenthion	0.05	19.1	0.0047	0.0125	13.9	90	278.0 → 109.0	15	278.0 → 169.0	15
17	Imazalil	1	24.1	0.0607	0.2500	9.4	86	214.9 → 173.0	5	216.8 → 175.0 172.9 → 145.0	5 15

No.	Pesticide	MRL	RT (min)	LOD	LOQ	RSD (%)	REC (%)	Quant	CE (V)	Qual	CE (V)
18	Kresoxim-methyl	0.2	25.0	0.0184	0.0500	13.9	88	116.0 → 89.0	15	131.0 → 89.0 116.0 → 63.0	30 30
19	Methidathion	0.05*	22.6	0.0056	0.0125	14.7	102	144.9 → 85.0	5	144.9 → 58.1 85.0 → 58.0	15 5
20	Prochloraz	0.2	31.9	0.0134	0.0500	10.6	85	180.0 → 138.0	10	310.0 → 70.0 308.0 → 70.0	15 15
21	Pyraclostrobin	0.1	34.7	0.0094	0.0250	12.2	102	132.0 → 77.1	20	132.0 → 104.0 164.0 → 132.1	15 10
22	Quintozene	0.3	13.4	0.0006	0.0025	8.6	95	294.9 → 236.8	15	294.9 → 142.9 236.8 → 143.0	45 30
23	Tebuconazole	0.1	27.8	0.0084	0.0250	10.3	108	250.0 → 125.0	20	125.0 → 99.0 125.0 → 89.0	20 15
24	Tetraconazole	0.1	19.9	0.0101	0.0250	12.6	106	336.0 → 217.9	20	336.0 → 203.8 170.9 → 136.0	30 10
25	Thiabendazole	2	21.8	0.0199	0.5000	1.5	87	201.0 → 174.0	15	201.9 → 175.0 173.9 → 65.0	15 30
26	Tiametoxam	0.05	20.7	0.0018	0.0125	4.2	112	212.0 → 139.0	15	247.0 → 182.0 247.0 → 212.0	15 5
27	Trifloxystrobin	0.05	27.3	0.0051	0.0125	14.4	95	116.0 → 89.0	15	172.0 → 145.1 116.0 → 63.0	15 30
IS	Triphenyl phosphate	–	28.0	–	–	–	–	325.0 → 169.0	20	326.0 → 233.0	10
28	Vinclozolin	0.05*	16.8	0.0044	0.0125	13.6	86	187.0 → 124.0	20	211.9 → 172.0 284.9 → 212.0	15 15

Note: \* Pesticide not permitted on this crop (NPC)

The recovery study was carried out to determine method accuracy and precision. For each blank matrix, three levels at 1 MRL, ½ MRL, and ¼ MRL were prepared. For the pesticides that do not have an established MRL, it was set as 0.01 mg/kg, with 0.05 mg/kg for the compounds that have MRLs set by the European Union, as shown on Table 1.

MRMs of the compounds were selected based on the Agilent Pesticide and Environmental Pollutants MRM Database (G9250AA) and Agilent Pesticide Analysis Reference Guide [5].

The LOQs for the pesticides were determined based on the recovery and RSD results, and defined as the selected lowest-validated spiked level meeting the requirement described in SANCO/12571/2013 [1], where recovery should be 70 to 120% and RSD below 20%. Five replicates were used. The LODs were calculated as three times the RSD of the spiked samples at their assigned LOQ levels.

### Real world sample analysis, quantitative method

Tommy Atkins mangoes were purchased at local groceries. Twenty lots of samples (from different stores) were analyzed. Three mangoes from each lot were extracted. The samples were chopped (including peel), blended, homogenized, and extracted soon after.

Using the quantitative method, it was possible to verify the presence of azoxystrobin, cyhalothrin, cypermethrin, difenoconazole, imazalil, prochloraz, pyraclostrobin, tebuconazole, tetraconazole, and thiabendazole below MRL. We also found chlorpyrifos, thiamethoxam, trifloxystrobin, and vinclozolin, none of which are permitted on mango. Two samples had concentrations of prochloraz above the MRL established by ANVISA, showing levels of 0.85 and 0.92 mg/kg (MRL = 0.2 mg/kg). The full list of pesticides is in Table 2.

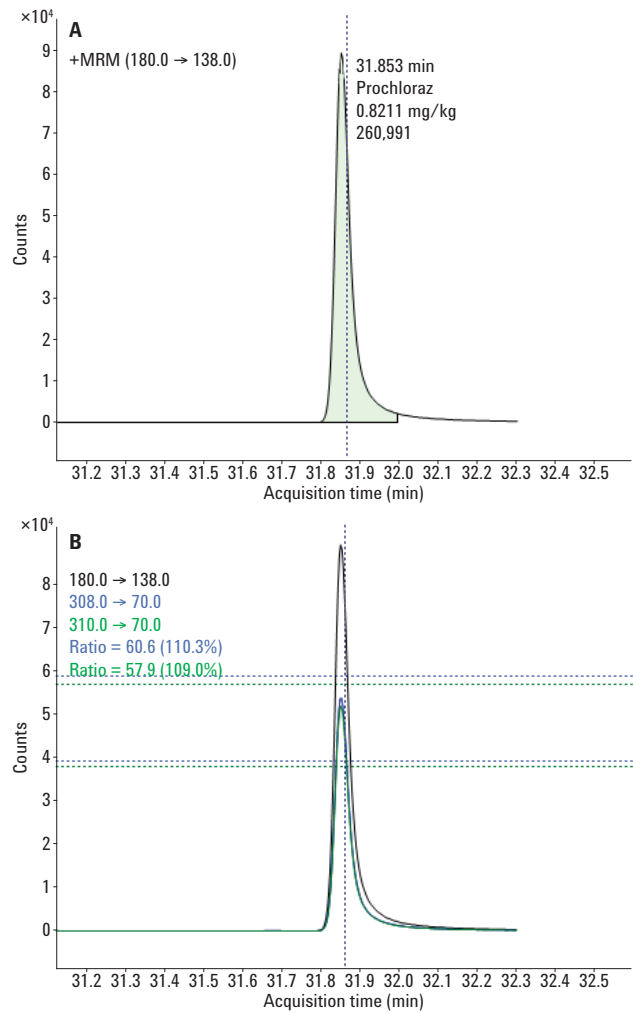


Figure 2. Prochloraz peak detected in a mango sample purchased at São Paulo supermarket.

Table 2. Pesticides detected in mango using the quantitative method.

Compound	Sample no.																			
	1	2	3	4	5	6	7	8	9	10	11	12	13	14	15	16	17	18	19	20
1 BHC- <i>alpha</i>	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-
2 Quintozene	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-
3 Diazinon	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-
4 Vinclozolin	-	-	-	-	-	-	-	<LOQ	<LOQ	-	<LOQ	<LOQ	-	-	-	-	-	-	-	-
5 Chlorpyrifos	-	-	-	-	-	-	-	<LOQ	-	-	-	-	-	-	-	<LOQ	-	-	-	-
6 Fenthion	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-
7 Tetraconazole	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	<LOQ	-
8 Tiametoxam	-	-	-	-	-	-	-	-	-	-	0.049	-	-	-	0.023	-	-	-	-	-
9 Thiabendazole	<LOQ	-	-	-	-	-	<LOQ	-	-	-	-	-	-	-	-	-	-	-	-	-
10 Methidathion	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-
11 Endosulfan I	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-
12 Imazalil	<LOQ	-	-	-	<LOQ	<LOQ	<LOQ	<LOQ	-	-	<LOQ	<LOQ	-	-	<LOQ	-	-	-	-	-
13 Kresoxim-methyl	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-
14 Endosulfan II	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-
15 Ethion	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-
16 Endosulfan sulfate	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-
17 Trifloxystrobin	-	-	-	-	-	-	-	-	-	-	-	<LOQ	-	<LOQ	<LOQ	-	-	-	-	-
18 Tebuconazole	-	-	0.0817	<LOQ	<LOQ	-	-	-	-	-	-	<LOQ	-	<LOQ	<LOQ	<LOQ	<LOQ	<LOQ	<LOQ	-
19 Carbosulfan	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-
20 Bifenthrin	-	-	-	-	-	-	-	-	-	-	-	-	-	-	<LOQ	<LOQ	-	-	-	-
21 Cyhalothrin ( $\lambda$ )	-	-	-	-	-	-	<LOQ	<LOQ	-	-	<LOQ	-	-	<LOQ	<LOQ	-	-	<LOQ	<LOQ	<LOQ
22 Prochloraz	-	-	-	-	-	-	<LOQ	0.92	-	-	-	-	-	-	-	0.8501	<LOQ	0.114	-	-
23 Cypermethrin I	-	-	-	<LOQ	-	-	-	-	-	-	-	<LOQ	-	<LOQ	<LOQ	-	-	-	<LOQ	-
24 Etofenprox	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-
25 Pyraclostrobin	-	0.0212	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	<LOQ	-	-
26 Difenconazole I	-	-	-	-	-	-	-	-	-	-	<LOQ	<LOQ	-	-	<LOQ	0.2225	<LOQ	0.0604	-	<LOQ
27 Azoxystrobin	-	-	<LOQ	<LOQ	<LOQ	-	-	-	-	-	-	-	-	<LOQ	<LOQ	<LOQ	<LOQ	<LOQ	-	<LOQ
28 Famoxadone	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-

Note: <LOQ = when peak was detected but the value was below LOQ, and above LOD. Values highlighted in red were above MRL.

### Real world sample analysis, qualitative screening

The samples were also analyzed by the qualitative method, with even more pesticides detected (Table 3).

The concentrations were calculated by injecting specific standards after detection.

Table 3. Pesticides found in mango by the qualitative method.

Compound	Sample no.																			
	1	2	3	4	5	6	7	8	9	10	11	12	13	14	15	16	17	18	19	20
1 Cyfluthrin	-	-	-	-	<0.03	<0.03	0.05	0.08	<0.03	<0.03	<0.03	<0.03	<0.03	<0.03	-	-	-	-	-	-
2 Dimethoate	-	-	-	-	-	-	-	0.6	-	-	-	-	-	-	-	-	-	-	-	-
3 Epoxiconazole	-	-	-	-	<0.05	-	-	-	-	-	<0.05	-	-	-	-	-	-	-	-	<0.05
4 Fenoxaprop-ethyl	-	-	-	-	-	-	<0.01	-	-	-	-	-	-	-	-	-	-	-	-	-
5 Mirex	-	-	-	-	-	-	-	-	-	-	<0.01	<0.01	<0.01	<0.01	-	<0.01	-	-	-	-
6 Omethoate	-	-	-	-	-	-	-	0.5	-	-	-	-	-	-	-	-	-	-	-	-
7 Permethrin	-	-	-	-	-	-	-	-	<0.03	<0.03	<0.03	<0.03	<0.03	<0.03	-	-	-	-	-	-
8 Propoxur	-	-	-	-	-	-	-	0.01	-	-	-	-	-	-	-	-	-	-	-	-
9 Trichlorophenol 2,4,6	-	-	-	-	-	-	<0.05	<0.05	-	-	-	-	-	-	-	-	<0.05	<0.05	<0.05	-

The comprehensive method allowed the detection of more than 250 pesticides at one run for screening purposes (Table 4). This method was easily created using the G9250AA Pesticide and Environmental Pollutants MRM Database, since the retention times of the compounds are already listed for the Pesticide Analyzer 411 configuration. In the past, the

preparation of the calibration curve for this number of compounds is very troublesome, and the daily checking of method integrity becomes much more involved. Screening first, and preparing calibration curves of identified compounds afterwards, makes the process faster, easier to perform, and far more cost-effective.

Table 4. Pesticides included in the MRM qualitative screening method.

1	3 OH Carbofuran	45	Cinerina I	90	Disulfoton
2	Acephate	46	Cinerina II	91	Disulfoton sulfone
3	Acetamiprid	47	Clomazone	92	Diuron
4	Acrinathrin	48	Cyanazine	93	Diuron metabolite (3,4-dichlorophenyl isocyanate)
5	Alachlor	49	Cycloate	94	DMSA (dichlofluanid metabolite)
6	Aldrin	50	Cyfluthrin I	95	DMST (tolylfluanid metabolite)
7	Ametryn	51	Cyfluthrin II (CAS no. 68359-37-5)	96	Endosulfan I ( <i>alpha</i> isomer)
8	Amitraz	52	Cyfluthrin III (CAS no. 68359-37-5)	97	Endosulfan II ( <i>beta</i> isomer)
9	Anilazine	53	Cyfluthrin IV (CAS no. 68359-37-5)	98	Endosulfan sulfate
10	Atrazine	54	Cyhalothrin ( <i>gamma</i> )	99	Endrin
11	Atrazine-desethyl	55	Cyhalothrin ( <i>lambda</i> )	100	Endrin aldehyde
12	Azinphos-ethyl	56	Cypermethrin I	101	Endrin ketone
13	Azinphos-methyl	57	Cypermethrin II	102	Epoxiconazole
14	Azoxystrobin	58	Cypermethrin III	103	EPTC
15	Benalaxyl	59	Cypermethrin IV	104	Ethion
16	Benfluralin	60	Cyproconazole	105	Ethylan (ethyl-DDD, Perthane)
17	Benfuracarb	61	Cyromazine	106	Etofenprox (ethofenprox)
18	BHC- <i>alpha</i> (benzene hexachloride)	62	DDD- <i>o,p'</i>	107	Famoxadone
19	BHC- <i>beta</i>	63	DDD- <i>p,p'</i>	108	Fenamidone
20	BHC- <i>delta</i>	64	DDE- <i>o,p'</i>	109	Fenamiphos (phenamiphos)
21	BHC- <i>epsilon</i>	65	DDE- <i>p,p'</i>	110	Fenamiphos sulfone
22	BHC- <i>gamma</i> (lindane, <i>gamma</i> HCH)	66	DDT- <i>o,p'</i>	111	Fenarimol
23	Bifenthrin	67	DDT- <i>p,p'</i>	112	Fenchlorphos oxon
24	Bromopropylate	68	Deltamethrin	113	Fenitrothion
25	Butylate	69	Demeton-S	114	Fenoprop-methyl
26	Captafol	70	Demeton-S-methyl	115	Fenoxaprop-ethyl
27	Captan	71	Demeton-S-methyl sulfone	116	Fenoxaprop- <i>p</i> -ethyl
28	Carbaryl	72	Diazinon	117	Fenpropathrin
29	Carbofuran	73	Diazinon-oxon (diazoxon)	118	Fenson
30	Carbofuran, 3-keto-	74	Dichlofluanid	119	Fensulfothion
31	Carbofuran, 7-phenol-	75	Dichlorvos	120	Fensulfothion sulfon
32	Carbophenothion	76	Dicloran (dichloran)	121	Fenthion
33	Carbophenothion-methyl (methyl trithionate)	77	Dicofol, <i>o,p'</i> -	122	Fenthion oxon sulfone
34	Carbosulfan	78	Dicofol, <i>p,p'</i> -	123	Fenthion sulfone
35	Chlordane- <i>cis</i> ( <i>alpha</i> )	79	Dicrototos (dicrototos)	124	Fenthion sulfoxide
36	Chlordane- <i>oxy</i>	80	Dieldrin	125	Fenvalerate I
37	Chlordane- <i>trans</i> ( <i>gamma</i> )	81	Difenoconazole I	126	Fenvalerate II (CAS no. 51630-58-1)
38	Chlorfenson	82	Difenoconazole II (CAS no. 119446-68-3)	127	Fipronil
39	Chlorfenvinphos	83	Dimethenamid	128	Fipronil sulfide
40	Chlorobenzilate	84	Dimethenamid-P	129	Fipronil sulfone
41	Chlorothalonil	85	Dimethoate	130	Fluazinam
42	Chlorpropham	86	Dinocap I	131	Flufenoxuron
43	Chlorpyrifos	87	Dinocap II	132	Flumetralin
44	Chlorpyrifos-methyl	88	Dinocap III	133	Fluquinconazole
		89	Dinocap IV		



134	Fluthiacet-methyl	175	Paraoxon	215	Prothioconazole-desthio
135	Flutriafol	176	Paraoxon-methyl	216	Pyraclostrobin
136	Folpet	177	Parathion	217	Pyrethrin I
137	Fonofos	178	Parathion-methyl	218	Pyrethrin II
138	Formothion	179	Pebulate	219	Pyridaphenthion
139	Heptachlor	180	Pendimethalin (penoxaline)	220	Pyriproxyfen
140	Heptachlor endo-epoxide (isomer A)	181	Pentachloroaniline	221	Quintozene
141	Heptachlor exo-epoxide (isomer B)	182	Pentachlorobenzene	222	Quintozene metabolite (pentachlorophenyl)
142	Hexachlorobenzene	183	Pentachlorophenol	223	Ronnel (fenchlorphos)
143	Hexazinone	184	Permethrin I	224	Simazine
144	Imazalil	185	Permethrin II ( <i>trans</i> )	225	Spirodiclofen
145	Iprodione	186	Pethoxamid	226	Spiromesifen
146	Iprovalicarb I	187	Petoxamida	227	Tebuconazole
147	Iprovalicarb II (CAS no. 140923-17-7)	188	Phenthoate	228	Tecnazene (TCNB)
148	Jasmolina I	189	Phorate	229	Tefluthrin, <i>cis</i> -
149	Jasmolina II	190	Phorate oxon sulfone	230	Terbufos
150	Kresoxim-methyl	191	Phorate sulfone	231	Terbufos sulfone
151	Lactofen	192	Phorate sulfoxide	232	Terbutylazine
152	Malaoxon (metabolite of malathion)	193	Phoratoxon	233	Terbutylazine-desethyl
153	Malathion	194	Phosalone	234	Terbutryn
154	MCPA-butoxyethyl	195	Phosmet	235	Tetrachlorvinphos, E-isomer
155	MCPA-methyl	196	Phosmet oxon	236	Tetraconazole
156	MCPB-methyl	197	Piperonyl butoxide	237	Tetradifon
157	Metalaxyl	198	Pirimicarb	238	Tetrasul
158	Methamidophos	199	Pirimiphos-ethyl	239	Thiabendazole
159	Methidathion	200	Pirimiphos-methyl	240	Thiametoxan
160	Methiocarb	201	Prochloraz	241	Thiazopyr
161	Methiocarb sulfone	202	Procymidone	242	Thiobencarb (benthiocarb)
162	Methiocarb sulfoxide	203	Profenofos	243	Thiometon
163	Methoxychlor olefin	204	Profluralin	251	Trichlorophenol, 2,4,6-
164	Methoxychlor, <i>o,p'</i> -	205	Prometon	252	Trifloxystrobin
165	Methoxychlor, <i>p,p'</i> -	206	Prometryn	253	Trifluralin
166	Mevinphos	207	Pronamide (propyzamide)	254	Trinexapaque
167	Mirex	208	Propaquizafop	255	Triphenyl phosphate
168	Molinate	209	Propargite	256	Vamidothion
169	Monocrotophos	210	Propazine	257	Vernolate
170	Naled	211	Propham	258	Vinclozolin
171	Nitrofen	212	Propiconazole I		
172	Omethoate	213	Propiconazole II (CAS no. 60207-90-1)		
173	Oxadiazon	214	Propoxur		
174	Oxamyl				

## Conclusions

This study developed two methods for pesticide residue analysis. The first was used to quantify pesticides that have MRLs determined by Brazilian legislation, with good recovery and detection limits. The second method was used to verify the presence of 258 pesticides and metabolites simultaneously. By having two methods, it is possible to simplify the preparation of daily calibration curves for

accurate quantification of listed pesticides, and also check if any pesticide not allowed for use with a particular crop or commodity has been used. For routine pesticide analysis, it is important to check as many compounds as possible. Our study showed that several banned pesticides were still being used, as well as pesticides that should not be used on mango.

## References

1. Anon. *Guidance Document on Analytical Quality Control and Validation Procedures for Pesticide Residues Analysis in Food and Feed, SANCO/12571/2013, 19 Nov 2014*. European Commission, Health and Consumer Protection Directorate-General, Brussels, Belgium.
2. Nanakaan Saave. *Export Factsheet – Mangoes*. International Trade Centre, Geneva, Switzerland, **2011**.
3. Araújo, José. *Mercado e Comercialização da Manga*. Embrapa, Brasília - DF, Brazil, **2004**.  
<http://sistemasdeproducao.cnptia.embrapa.br/FontesHT/ML/Manga/CultivodaMangueira/mercado.htm>
4. ANVISA. Brazilian Health Surveillance Agency, Brasília - DF, Brazil.  
<http://portal.ANVISA.gov.br/wps/content/ANVISA+Portal/ANVISA/Inicio/Agrotoxicos+e+Toxicologia>
5. Anon. *GC/MS/MS Pesticide Residue Analysis; Pesticide Analysis Reference Guide*, Agilent Technologies, Inc. Publication number 5991-2389EN, **2013**.
6. Anon. *Foods of Plant Origin—Determination of Pesticide Residues Using GC-MS and/or LC-MS/MS Following Acetonitrile Extraction/Partitioning and Clean-up by Dispersive SPE (QuEChERS-method)*. EN 15662 Version 2008. European Standards, Pilsen, Czech Republic.
7. Anon. *Specification for Agricultural Pesticides in Agriculture*. Food and Agriculture Organization of the United Nations, Rome, Italy.  
<http://www.fao.org/agriculture/crops/thematic-sitemap/theme/pests/jmps/ps-new/en/>
8. Anastassiades, M.; Lehotay, S. J. Fast and Easy multiresidue method employing acetonitrile extraction/partitioning and “Dispersive solid-phase extraction” for the determination of pesticide residues in produce. *J. AOAC Int.* **2003**, *86*, 412-431.
9. Ciscato, C. H. P.; Gebara, A. B.; Monteiro, S. H. Pesticide residue monitoring of Brazilian fruit for export 2006-2007. *Food Additives and Contaminants: Part B* **2009**, *2*, 140-145.
10. Ciscato, C. H. P.; Gebara, A. B.; Barbosa, C. M.; Souza, G. S.; Hamada, N.; Silva, D. L. Pesticide Residues Evaluation in Pulp, Juice and Nectar of Fruits. *Environment and Pollution* **2012**, *1*, 39-43.
11. Anon. *Plano Nacional de Controle de Resíduos e Contaminantes em Produtos de Origem Vegetal*. Ministério da Agricultura Pecuária e Abastecimento (MAPA), Brasília - DF, Brazil.  
<http://www.agricultura.gov.br/vegetal/qualidade-seguranca-alimentos-bebidas/alimentos/residuos-e-contaminantes>.

## For More Information

These data represent typical results. For more information on our products and services, visit our Web site at [www.agilent.com/chem](http://www.agilent.com/chem).

[www.agilent.com/chem](http://www.agilent.com/chem)

Agilent shall not be liable for errors contained herein or for incidental or consequential damages in connection with the furnishing, performance, or use of this material.

Information, descriptions, and specifications in this publication are subject to change without notice.

© Agilent Technologies, Inc., 2015  
Printed in the USA  
July 2, 2015  
5991-6054EN



**Agilent Technologies**