

# Meeting the European Commission Performance Criteria for the Use of Triple Quadrupole GC-MS/MS as a Confirmatory Method for PCDD/Fs in Food and Feed Samples

Cristian Cojocariu<sup>2</sup>, Esteban Abad Holgado<sup>1</sup>, Manuela Abalos<sup>2</sup> and Paul Silcock<sup>2</sup>.

<sup>1</sup>Spanish Council for Scientific Research (CSIC), Institute of Environmental Assessment and Water Research, Barcelona, Spain.

<sup>2</sup>Thermo Fisher Scientific, Runcorn, Cheshire, United Kingdom.

## OVERVIEW

### Purpose

In this work, the performance of the Thermo Scientific™ TSQ 8000™ Evo Triple Quadrupole GC-MS/MS system for the analysis of Polychlorinated dibenzo-p-dioxins (PCDDs) and Polychlorinated dibenzo-p-furans (PCDFs) was assessed. For this, both solvent standards and food and feed samples were used to evaluate the instrument performance against the criteria for PCDD/Fs confirmation. Additionally, a direct comparison of the results obtained from food and feed sample extracts using the TSQ 8000 Evo GC-MS/MS system with those from a GC-HRMS was made.

### Methods

PCDD/Fs were analyzed in the standards and matrix samples using a TSQ 8000 Evo Triple Quadrupole GC-MS/MS instrument coupled with a Thermo Scientific™ TRACE™ 1310 GC system. Sample introduction was performed with a Thermo Scientific™ TriPlus™ RSH autosampler, and compound separation was achieved on a Thermo Scientific™ TraceGOLD TG-5iMS 60 m x 0.25 mm I.D. x 0.25 µm film capillary column.

### Results

The selectivity, sensitivity and repeatability of the method developed was shown to be more than adequate; meeting or exceeding the requirements stipulated in the legislation.

## INTRODUCTION

Until recently, legislation in the European Union required the confirmation and quantification of dioxins PCDD/Fs in contaminated samples by gas chromatography/high resolution mass spectrometry (GC-HRMS) instruments, which is still considered the "gold standard" approach. This technique has the sensitivity and specificity to be used for low level background monitoring as well as maximum and action levels in food and feed. However, recent advances in gas chromatography/triple-quadrupole mass spectrometry (GC-MS/MS) technology have allowed high sensitivity and selectivity to be achieved. These improvements have led to GC-MS/MS being considered a reliable tool that can be used to control the maximum levels for PCDD/Fs in food and feed as a full confirmatory method.<sup>4</sup> According to the revised EU regulation, when using GC-MS/MS, the following specific performance criteria for confirmation with GC-MS/MS technology should be fulfilled in addition to the criteria described previously by the European Commission<sup>1,3</sup>, except the obligation to use GC-HRMS:

- Resolution for each quadrupole to be set equal to or better than unit mass resolution (unit mass resolution defined as sufficient resolution to separate two peaks with one mass unit apart).
- Two specific precursor ions, each with one specific corresponding transition product ion for all labelled and unlabelled analytes should be used.
- Maximum permitted tolerance of relative ion intensities of  $\pm 15\%$  for selected transitions in comparison to calculated or measured values (average from calibration standards), applying identical MS/MS conditions, in particular collision energy and collision gas pressure, for each transition of an analyte.

## METHOD

### Sample Preparation

PCDD/F standards (EPA Method 1613 1613CSL to 1613CS4) containing the native and the <sup>13</sup>C-labelled compounds were obtained from Wellington Laboratories Inc. The following food and feed extracted samples were provided by the Institute of Environmental Assessment and Water Research, CSIC Barcelona, Spain 3x dry fish samples (previously used in inter-laboratory studies), one feed sample (internal reference material), and one milk powder sample (certified reference material).

Extraction and clean-up of the matrix samples was performed either by PowerPrep™ SPE system (feed sample) or by using a manual clean-up with multi-layer silica, followed by basic alumina and a final carbon column (milk, and fish samples). Final extracts were prepared in nonane.

The GC method setup is described in Table 1 below, while the MS method setup is described in Tables 2 and 3.

Tables 1 & 2. Instrument setup.

Trace 1310 GC Parameters		TSQ 8000 Evo Mass Spectrometer Parameters	
Injection Volume (µL):	2	Transfer line (°C):	280
Liner:	S51 Single Taper 4mm ID 78.5mm	Ionization type:	EI
Inlet (°C):	260	Ion source (°C):	300
Inlet Module and Mode:	Splitless	Electron energy (eV):	40
Carrier Gas, (mL/min):	He, 1.2	Acquisition Mode:	SRM
Oven Temperature Program:		O2 Gas Pressure (argon)(psi):	60
Temperature 1 (°C):	100	Collision Energy (eV):	see Table 3
Hold Time (min):	2	Q1 Peak Width (AMU):	0.7
Temperature 2 (°C):	250	Q3 Peak Width (AMU):	0.7
Rate (°C/min):	25		
Hold Time (min):	2		
Temperature 3 (°C):	285		
Rate (°C/min):	2.5		
Temperature 4 (°C):	330		
Rate (°C/min):	10		
Hold Time (min):	5		

Table 3. SRM transitions used for native and <sup>13</sup>C-labelled PCDD/Fs.

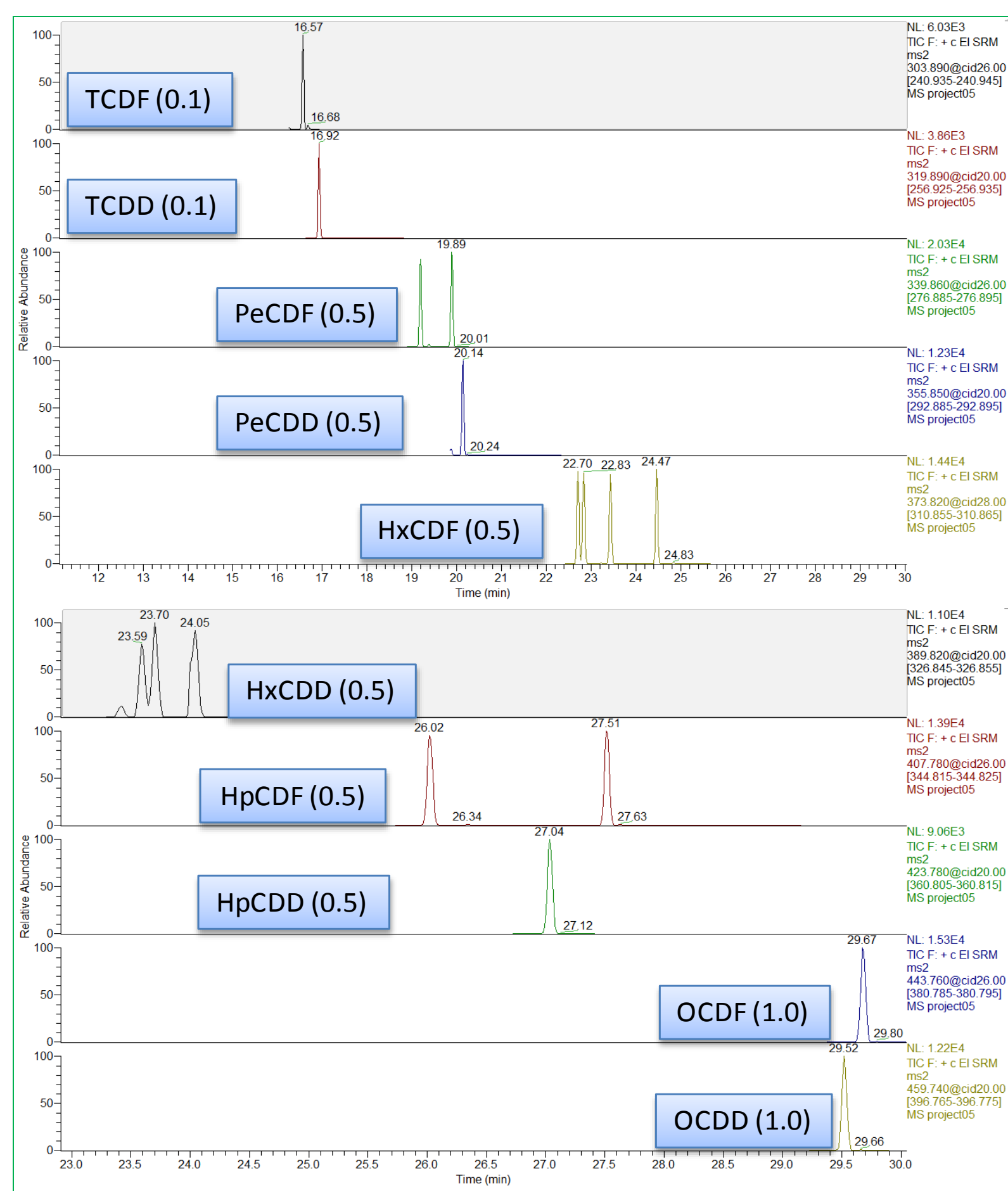
SRM Compound Information				SRM Compound Information			
Compound Name	Quant	Precursor Ion [Da]	Product Ion [Da]	Compound Name	Quant	Precursor Ion [Da]	Product Ion [Da]
<sup>13</sup> C-TCDF	Quant	315.94	251.97	<sup>13</sup> C-HxCDF	Quant	399.86	335.89
<sup>13</sup> C-TCDF	Quant	317.94	253.97	<sup>13</sup> C-HxCDF	Quant	401.86	337.89
TCDF	Quant	303.89	240.94	HxCDF	Quant	387.82	324.86
TCDF	Quant	305.89	242.94	HxCDF	Quant	389.82	326.86
<sup>13</sup> C-TCDD	Quant	331.94	267.97	<sup>13</sup> C-HpCDF	Quant	419.82	355.86
<sup>13</sup> C-TCDD	Quant	333.94	269.97	<sup>13</sup> C-HpCDF	Quant	421.82	357.86
TCDD	Quant	319.89	256.93	HpCDF	Quant	407.78	344.82
TCDD	Quant	321.89	258.93	HpCDF	Quant	409.78	346.82
<sup>13</sup> C-PeCDF	Quant	351.89	287.93	<sup>13</sup> C-HpCDD	Quant	435.82	371.85
<sup>13</sup> C-PeCDF	Quant	353.89	289.93	<sup>13</sup> C-HpCDD	Quant	437.81	373.85
PeCDF	Quant	339.86	276.89	HpCDD	Quant	423.78	360.81
PeCDF	Quant	341.86	278.89	HpCDD	Quant	425.77	362.81
<sup>13</sup> C-PCDD	Quant	367.89	303.93	<sup>13</sup> C-OCDD	Quant	469.78	405.81
<sup>13</sup> C-PCDD	Quant	369.89	305.93	<sup>13</sup> C-OCDD	Quant	471.78	407.81
PCDD	Quant	355.85	292.89	OCDD	Quant	457.74	394.77
PCDD	Quant	357.85	294.89	OCDD	Quant	459.74	396.77
<sup>13</sup> C-HxCDF	Quant	383.86	319.89	OCDF	Quant	441.76	378.79
<sup>13</sup> C-HxCDF	Quant	385.86	321.89	OCDF	Quant	443.76	380.79
HxCDF	Quant	371.82	308.86				
HxCDF	Quant	373.82	310.86				

## RESULTS

### Chromatography

Achieving sufficient chromatographic separation of the PCDD/F isomers is critical (sufficient defined as <25% peak to peak between 123478-HxCDF and 123678-HxCDF)<sup>1</sup>. Chromatography of PCDD/Fs was assessed in the lowest calibration standard (EPA1613-CLS) containing 0.1 pg/µL TCDD/F, 0.5 pg/µL PeCDD/F - HpCDD/F and 1.0 pg/µL OCDD/F. All the native and their corresponding <sup>13</sup>C-labelled internal standards were easily detected, excellent peak shape was obtained for all compounds (Figure 1), and 5% valley separation was achieved for HxCDF isomers (Figure 2).

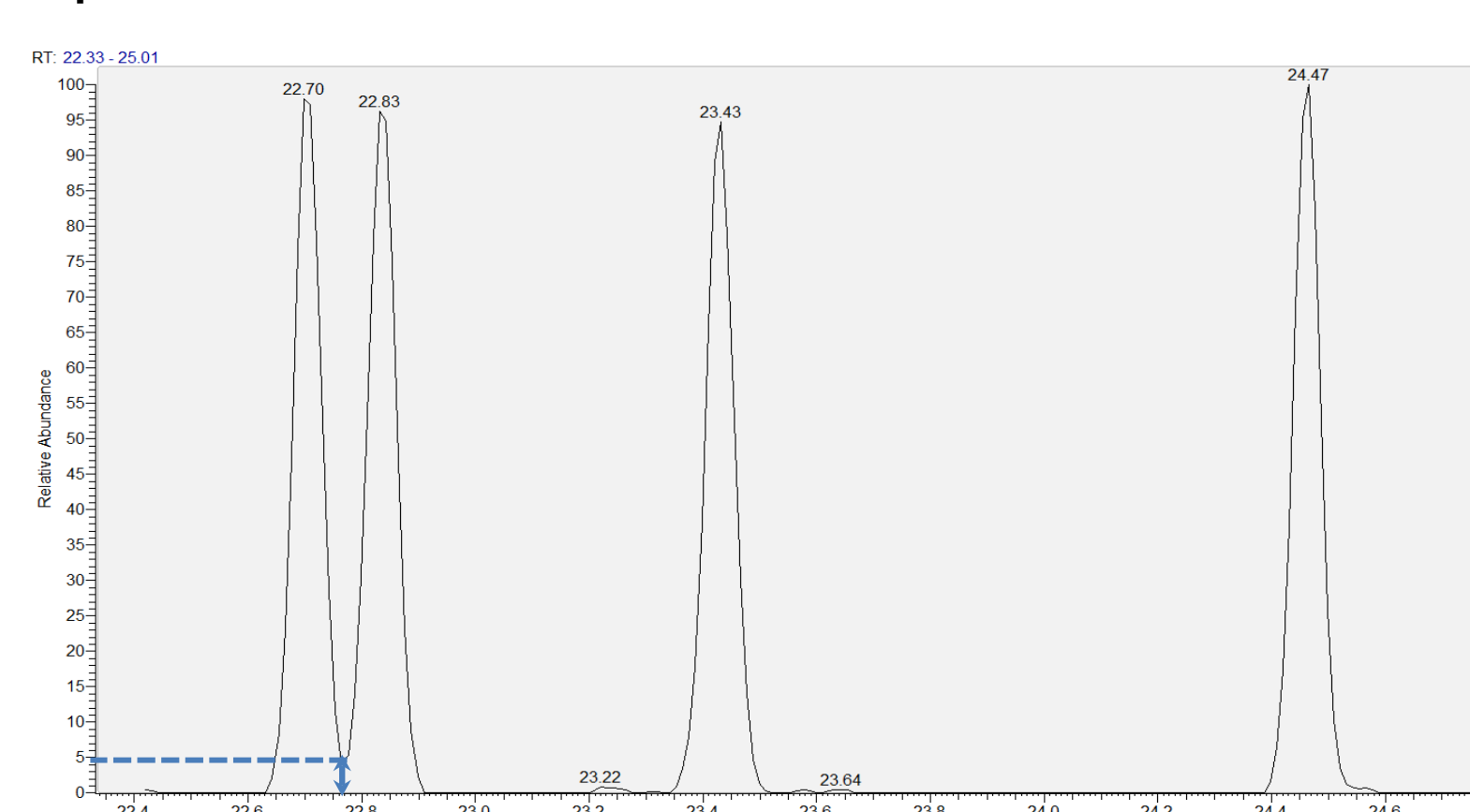
Figure 1. Chromatographic separation of native PCDD/Fs in the lowest standard (in brackets concentration in pg/µL). One SRM transition (quantification ion) per compound is shown.



### Data Analysis

Data processing was performed using the Thermo Scientific™ TargetQuan 3 software, which is specifically designed for the analysis of POPs using isotope dilution. The software streamlines quantitation based upon relative response factors (or optionally average responses), incorporates toxic equivalence factors (TEFs) to automatically calculate toxic equivalence quotients (TEQs) and finally total TEQ.

Figure 2. GC separation of HxCDF isomers showing 5% valley separation between 123478-HxCDF and 123678-HxCDF.

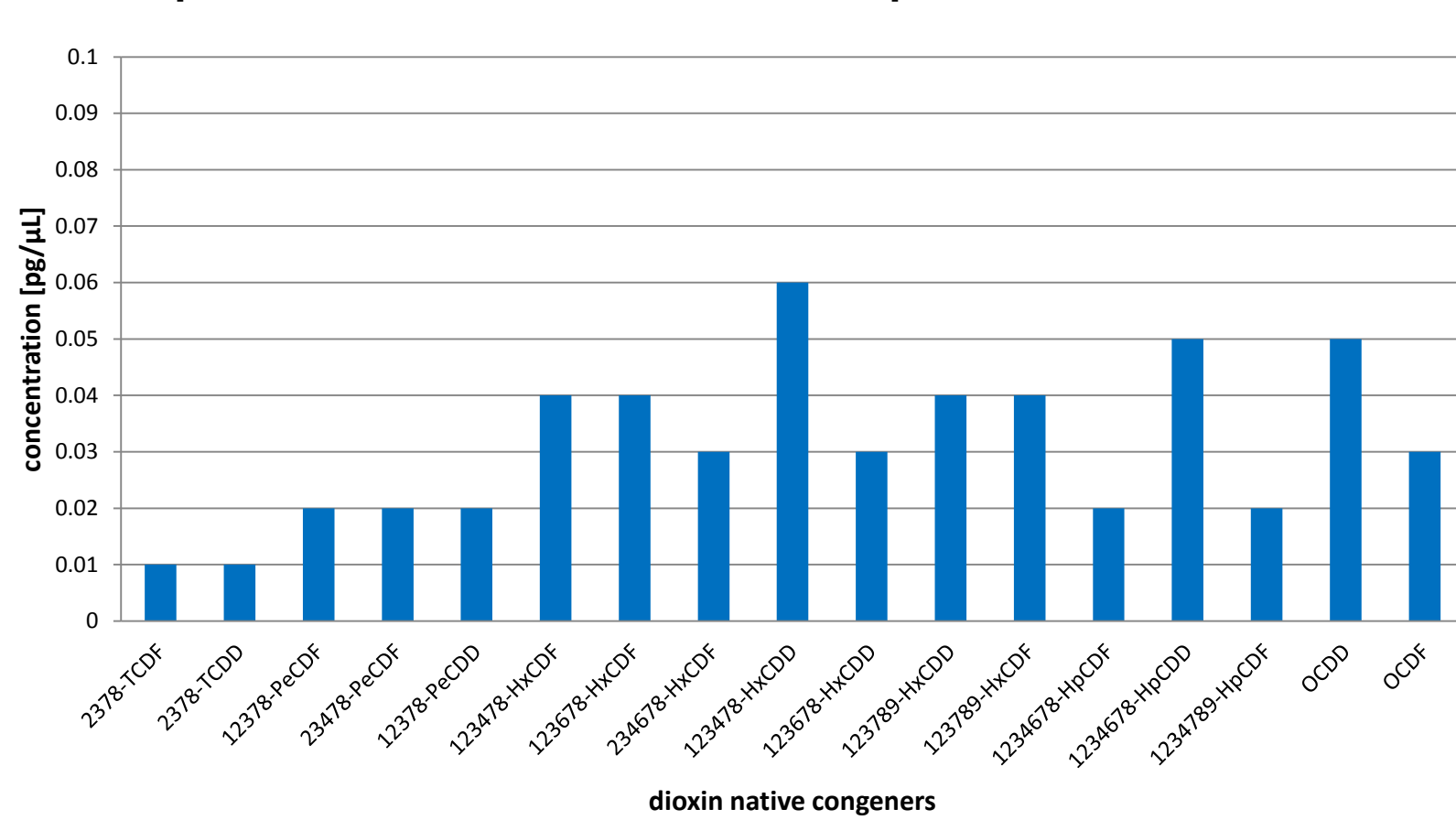


### Reaching the level of interest

For the analysis of PCDD/Fs, reaching the required limits of detection is critical to implementing a routine monitoring method. The limit of quantification (LOQ) for a confirmatory method should be about one fifth of the maximum level.<sup>1,3,5</sup>

The instrument LOQ was assessed by repeatedly (n = 10) injecting the lowest calibration standard (1613CSL) and three subsequent serial dilutions of this standard. The results of this test show that the LOQs for the PCDD/Fs analyzed were between 0.01 – 0.06 pg/µL (ion ratios and response factors RF at these levels still within  $\pm 15\%$  limit, % recovery of <sup>13</sup>C-labelled within the 60 – 120% limit) (Figure 3). The results of this experiment demonstrate that the TSQ 8000 Evo GC-MS/MS system can detect and confirm PCDD/Fs at low femtogram levels, thus meeting the detection limit requirements.<sup>3</sup>

Figure 3. LOQ calculation for PCDD/Fs from repeat injections of a serial dilution. Data indicate the LOQ for each congener with ion ratios and response factors values within the expected limits.



### Linearity of response

PCDD/F quantification is based on isotope dilution and uses RF type calibration where the average response factor of all the standards from an external calibration curve are taken into account to quantify the 17 toxic congeners.<sup>1,6</sup> Average RF %RSD values were calculated from duplicate measurements of a six point calibration curve measured at the beginning and at the end of the sample batch. The results of this experiment show excellent %RSD for all measured compounds with values between 0.8 – 3.6 %, well within the 15% limits established by EPA<sup>6</sup> (Table 4).

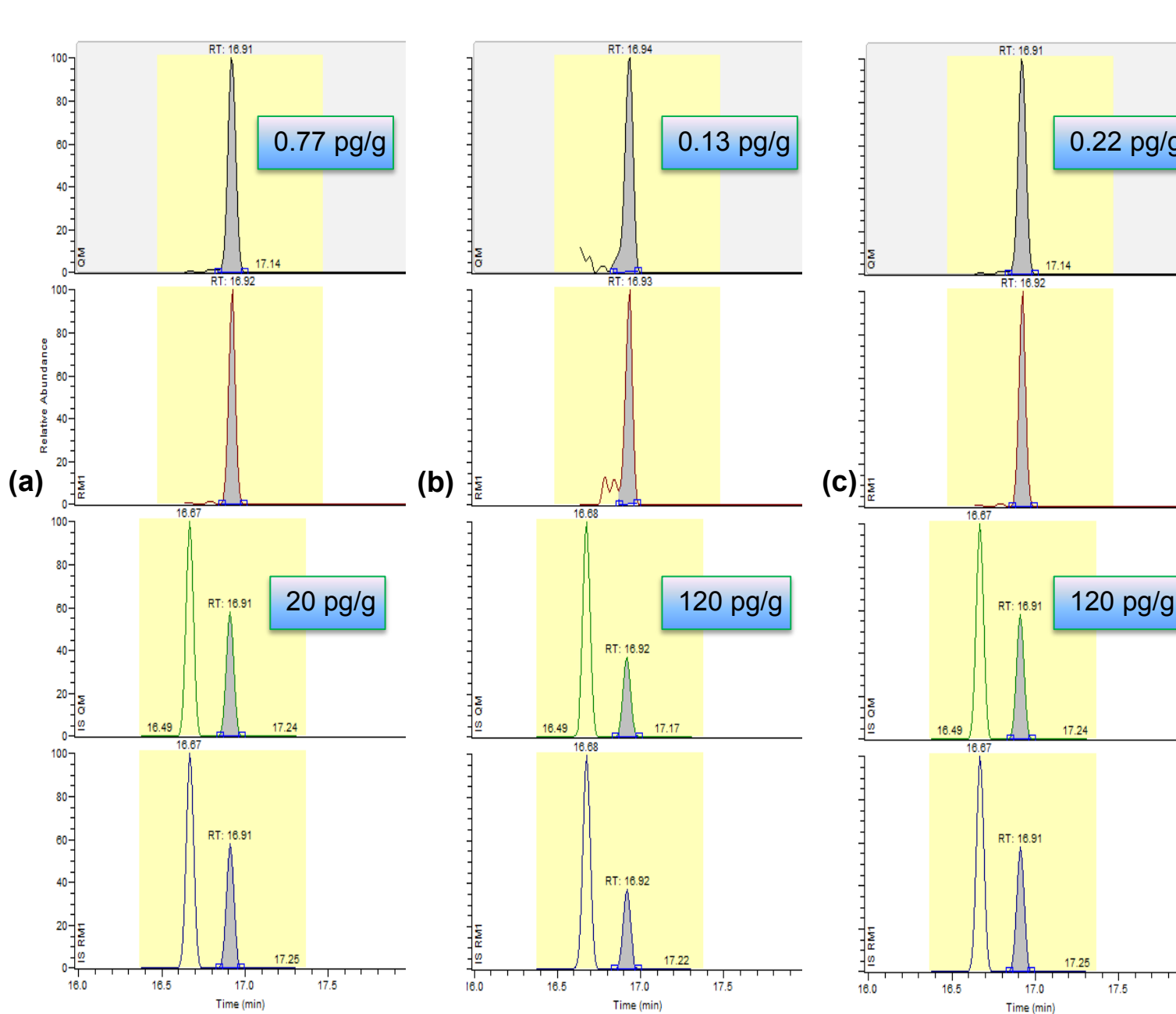
Table 4. Linearity of PCDD/Fs across six point calibration curve. The precision on the average response factor (%RSD) for each native compound is shown. Values represent duplicate measurements of each calibration point, measured at the beginning and end of a batch.

Compound	Concentration range (pg/µL)	Linearity / Calibration		
		Average RF	stdev	RF % RSD
2378-TCDF	0.1 - 40	1.04	0.02	1.9
2378-TCDD	0.1 - 40	1.12	0.02	2.2
12378-PeCDF	0.5 - 200	1.01	0.02	1.5
23478-PeCDF	0.5 - 200	1.03	0.02	1.6
12378-PeCDD	0.5 - 200	1.08	0.01	1.4
123478-HxCDF	0.5 - 200	1.03	0.01	1.2
123678-HxCDF	0.5 - 200	1.02	0.01	1.4
234678-HxCDF	0.5 - 200	1.06	0.03	3.2
123478-HxCDD	0.5 - 200	0.96	0.01	1.6
123678-HxCDD	0.5 - 200	1.21	0.04	3.6
123789-HxCDD	0.5 - 200	1.21	0.04	3.6
123789-HxCDF	0.5 - 200	1.36	0.11	0.8
1234678-HpCDF	0.5 - 200	1.06	0.02	1.7
1234678-HpCDD	0.5 - 200	1.07	0.02	2.1
1234789-HpCDF	0.5 - 200	1.12	0.02	2.2
OCDD	1.0 - 400	1.54	0.04	2.4
OCDF	1.0 - 400	1.09	0.03	3.2

### Quantification of PCDD/Fs in sample extracts

PCDD/Fs were quantified in the sample extracts prepared at CSIC, Barcelona. Excellent chromatographic separation with little matrix interference was observed for all native congeners for all sample analyzed. An example of chromatography is shown below for 2378-TCDD (Figure 4).

Figure 4. Example of chromatographic separation of 2,3,7,8-TCDD and its internal standard <sup>13</sup>C-2,3,7,8-TCDD present in the fish (a) feed (b) milk powder (c) samples. Calculated concentration (pg/g) is indicated.



PCDD/F content of each sample, expressed as WHO-TEQ pg/g, was determined for each sample analyzed and the results were compared with the existing data obtained for the same samples from the GC-HRMS. The calculated concentrations of each individual PCDD/F congener (as TEQ pg/g) were compared with the values obtained from the GC-HRMS. The data shows excellent agreement between the results obtained using the TSQ 8000 Evo GC-MS/MS system and that obtained using GC-HRMS (Figures 5 and 6). The concentrations of the PCDD/Fs in the samples cover a wide range. But importantly, there is always good agreement with the GC-HRMS results.

Figure 5. Individual contribution of each PCDD/F congener to the feed sample dioxin content (as TEQ pg/g) and comparison of TSQ 8000 Evo GC-MS/MS system results with the GC-HRMS values.

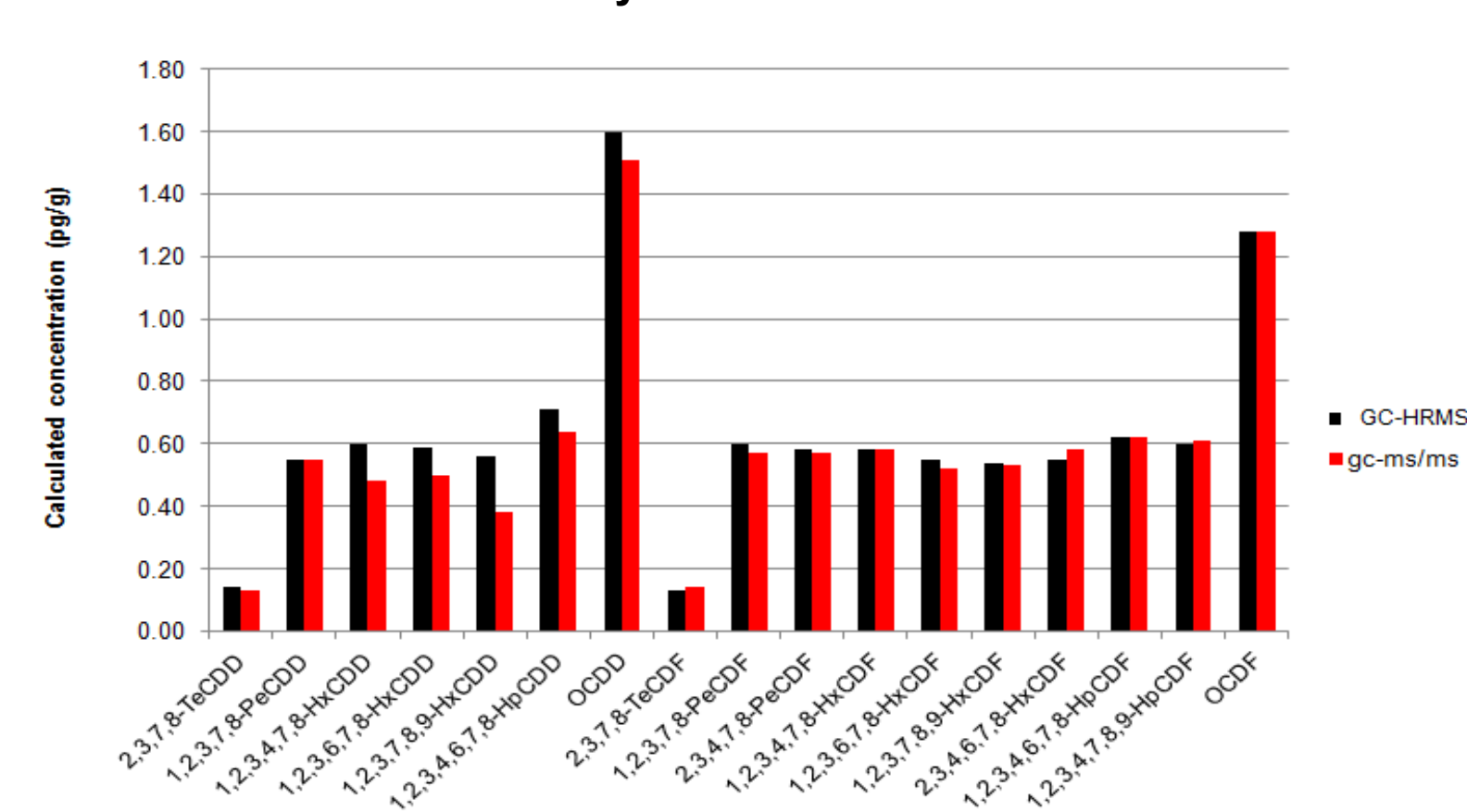
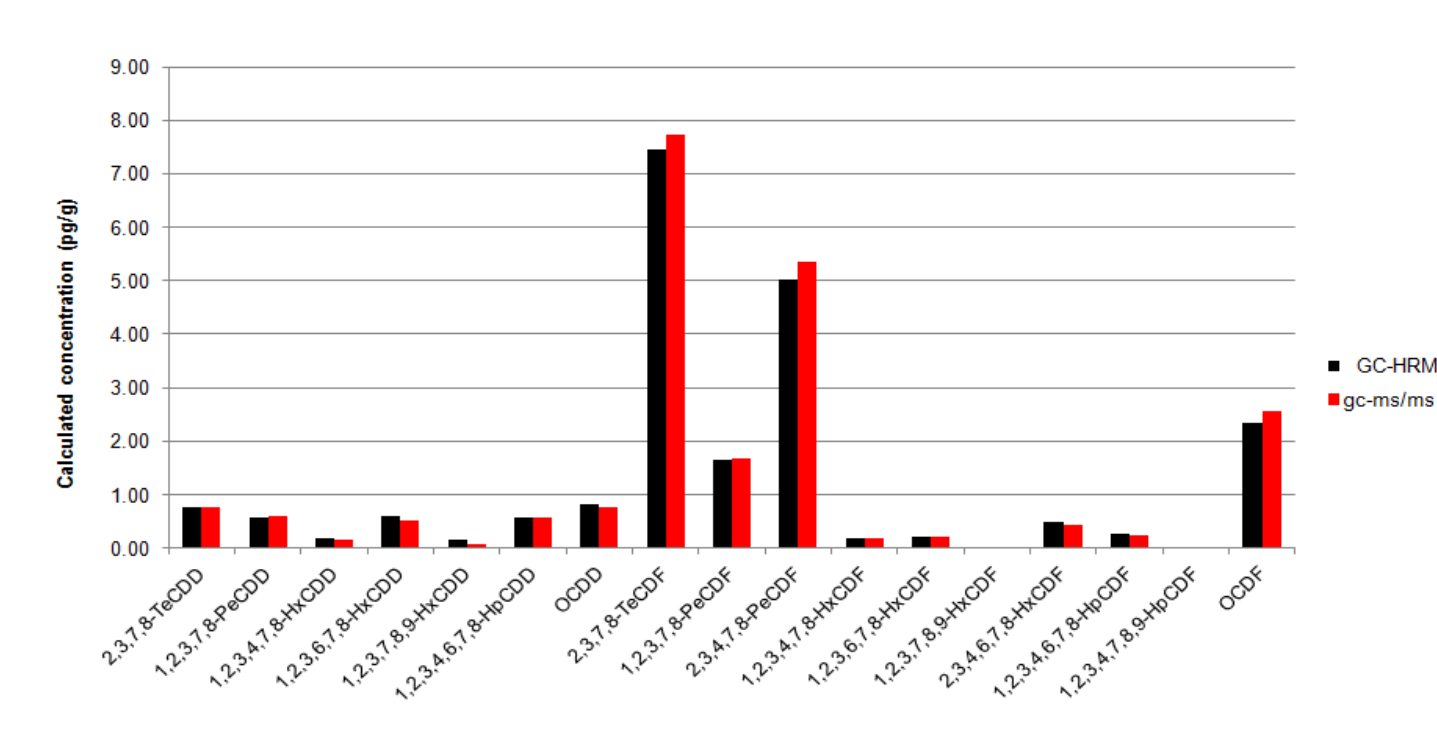


Figure 6. Individual contribution of each PCDD/F congener to the feed sample dioxin content (as TEQ pg/g) and comparison of the TSQ 8000 Evo GC-MS/MS system results with the GC-HRMS values.

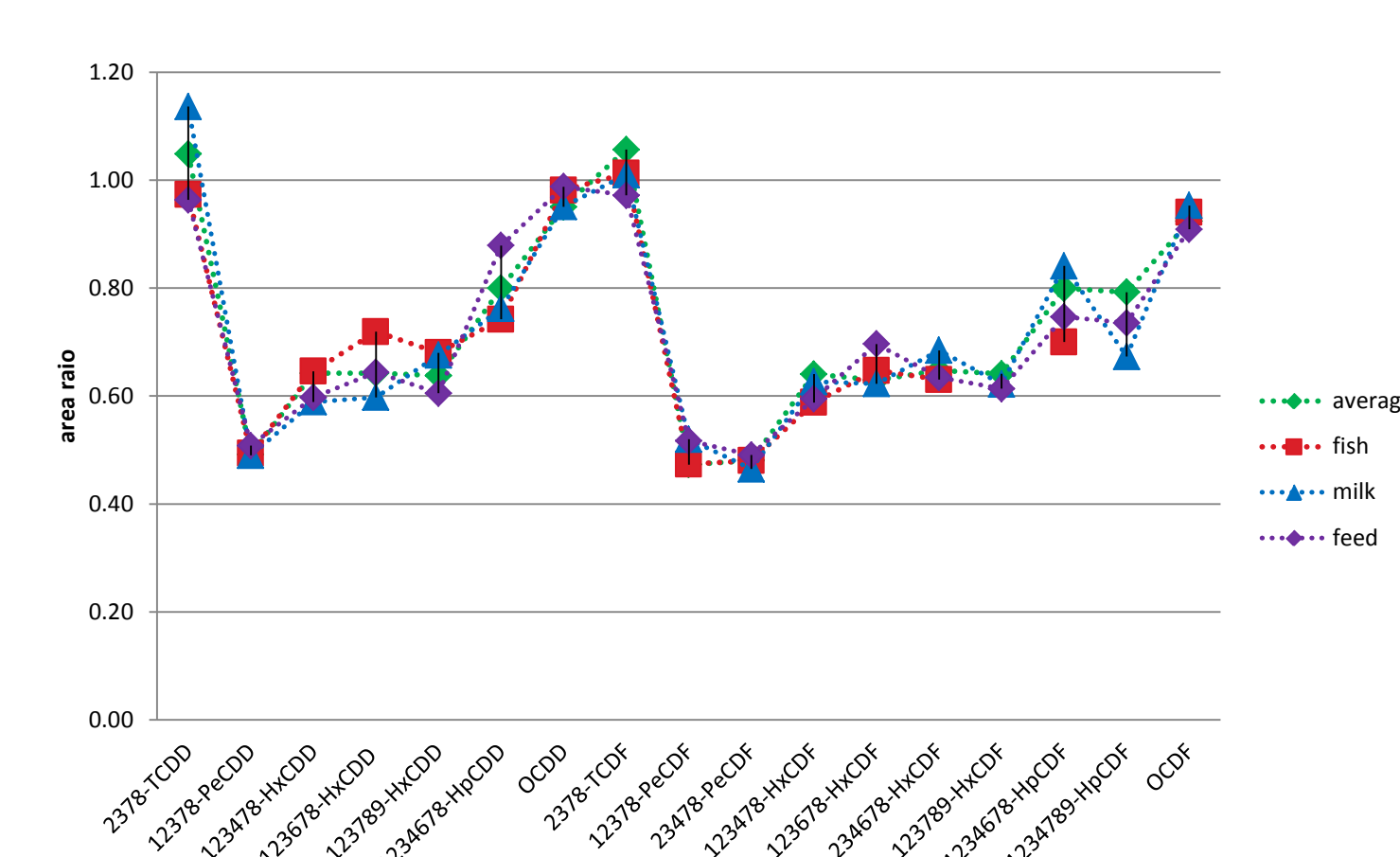


The total PCDD/F content of each sample obtained from the TSQ 8000 Evo GC-MS/MS system analysis was plotted against the sector instrument data, with the calculated deviation not exceeding 5%.

### Ion ratio abundance

The ion ratio (IR) abundance for selected transitions of each of the 17 PCDD/F congeners was measured in each of the samples analyzed and the values compared with the measured ion ratio values (average from calibration standards CSL-CS4). The results of this experiment show that all the IR for the compounds analyzed were within the 15% tolerance meeting the EU criteria for dioxin confirmation<sup>2</sup> (Figure 7).

Figure 7. Comparison of the ion ratio abundance of each of the 17 PCDD/F in the samples extracts with the average IR values derived from the calibration standards (CSL-CS4).



## CONCLUSIONS

Taken together, the results of this evaluation demonstrates that the TSQ 8000 Evo GC-MS/MS system is an extremely effective tool for routine analysis of PCDD/Fs meeting all the European Commission requirements for the confirmation of dioxins in food and feed samples. The results obtained with the TSQ 8000 Evo GC-MS/MS instrument demonstrate that this is a highly sensitive and selective analytical system that can be confidently used for PCDD/Fs detection and confirmation in food and feed samples.

- The TSQ 8000 Evo GC-MS/MS system, together with the TRACE 1310 GC system and the TargetQuan 3 data processing and reporting software, constitute a comprehensive system solution for dioxin and furan analysis in complex samples.
- Excellent reproducibility, linearity, sensitivity, and selectivity were obtained in all the experiments performed with standards and sample extracts.
- Moreover, the calculated PCDD/Fs TEQ values for the matrix samples were in very good agreement with those derived from the sector instrument, the results recommending this system for routine and confident analysis of PCDD/Fs.

## REFERENCES

- European Commission, Commission Regulation No 1883, Off. J. Eur. Union, L 364, 32-43, 2006.
- European Commission, Commission Regulation No 1259, Off. J. Eur. Union, L 320, 18-23, 2011.
- European Commission, Commission Regulation No 252, Off. J. Eur. Union, L 84, 1-22, 2012.
- European Commission, Commission Regulation No 589, Off. J. Eur. Union, L164, 18-40, 2014.
- Kotz A., Malisch R., Focant J., Eppe G., Cederberg Tommy Licht, Rantakokko P., Fürst P., Bernsmann T., Leondiadis L., Lovasz C., Scortichini G., Dielti G., di Domenico A., Ingelido A. M., Traag W., Smith F., Fernandes A., Organohalogen Compounds, Vol. 74, 156-159, 2012.
- US EPA Method 1613: Tetra-through octa-chlorinated dioxins and furans by isotope dilution HRGC/HRMS (Revision B), 1994.

## TRADEMARKS/LICENSING

© 2016 Thermo Fisher Scientific Inc. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. PowerPrep is a trademark of FMS, Inc. This information is not intended to encourage use of these products in any manner that might infringe the intellectual property rights of others.

**ThermoFisher**  
SCIENTIFIC