

Enhancement of Linearity and Response in Charged Aerosol Detection

Christopher Crafts, Marc Plante, Bruce Bailey, Ian Acworth, Thermo Fisher Scientific, Chelmsford, MA, USA



Overview

Purpose: This work evaluates the application of a power transformation algorithm during data acquisition with a charged aerosol detection. Its effect on increased linear range, peak shape, response, resolution, signal-to-noise and limits of detection will be discussed.

Methods: The effect of changing the power function variable was examined using standards ranging from nanogram to mid-microgram amounts on column. It was also evaluated using a range of analytes showing different physicochemical properties and chromatographic chemistries.

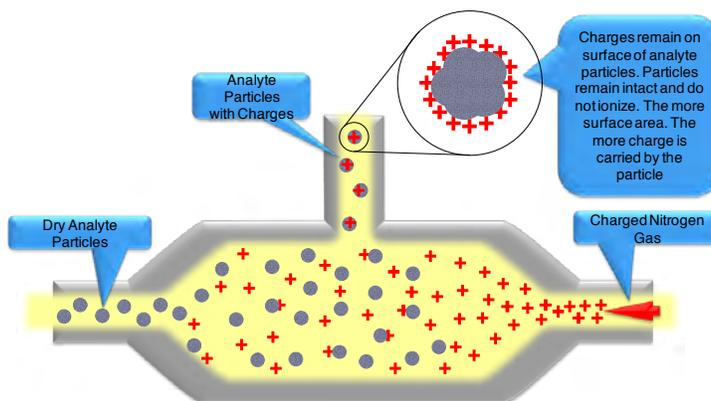
Results: Linearity was improved over the mass range analyzed with correlation coefficients generally >0.999 or greater. Significant improvements in signal-to-noise and resolution along with decreases in peak asymmetry and peak width were also observed but were dependent upon the algorithm value used.

Introduction

Raw data collected by analytical detectors rarely correlates directly to the data points that are analyzed by the chromatographic software. Rather the data has already been manipulated in some way within the detector, typically post data collection. The most common examples are filtering and smoothing algorithms. The power transformation is simply another internal data manipulation aimed at improving the outputted data quality without changing the true separation. The use of power transformation can often improve the appearance of non-linear detectors. While these settings are typically not fully disclosed or described to the end user, they may be implemented to improve the appearance or linearity of the results.

The Thermo Scientific Dionex Corona ultra RS Charged Aerosol Detector (CAD™) is a highly sensitive universal detector that can deliver near uniform response for non-volatile analytes.¹ Charged aerosol detection works by nebulizing the HPLC eluent into fine droplets that are then dried to form particles. The size of these particles directly correlates to the mass injected. After drying particles then collide with a stream of charged nitrogen and the charge is transferred to the surface of the particle (Figure 1). The charge on the particles continue is collected and converted to current using a highly sensitive electrometer. Charged aerosol detection is non-linear as the amount of analyte being measured depends on the charge that can be placed on the surface of the particle. This in turn is dependent upon the volume of the particle being formed. The relationship between particle surface area and volume is not linear. This is observed experimentally as the response curves for the Corona™ ultra RS™ Charged Aerosol Detector are typically expressed as the 2nd order polynomial function over a range of up to 4 orders of magnitude.² The application of power function can now correct for this physical characteristic and deliver a more linear response over a larger dynamic range.

FIGURE 1: Mixing Chamber of the Corona ultra RS Charged Aerosol Detector.



A power function is a simple mathematic transformation where the individual data points are raised to a preset value. The possible values are between 1.0 and 2.0 for the Corona ultra RS detector. This power function value (PFV) can vary depending on the chemistry in use and the mass range being analyzed. While its primary function is to increase linear range, several other benefits may result from a power transformation^{2,3,4}. Data are presented to demonstrate the impact of the power function on linearity, peak width, resolution, and signal-to-noise (S/N).

Methods

Sample Preparation

Standards of 14 compounds were dissolved in appropriate solvents at concentration of ~ 2 mg/mL. These compounds were then diluted using serial dilution. The mass range tested was from ~50 ng to as high as 14 μg on-column (7 to 10 point response curves) depending on the analysis in question. Further detail on the analysis range can be found in the figures below.

Liquid Chromatography

Thermo Scientific Dionex UltiMate 3000 RSLC system along with a Corona ultra RS Charged Aerosol Detector was used for all data generation.

Exact analytical conditions varied for the different compounds analyzed and are not relevant to the data presented here. Method conditions are available upon request.

Data Analysis

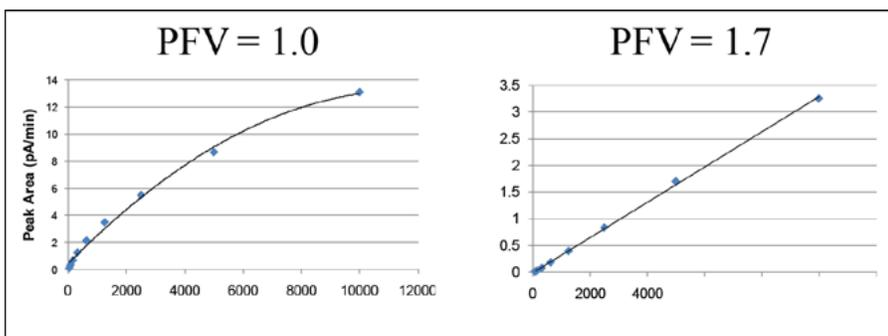
Thermo Scientific Chromeleon Chromatography Data System 6.8 SR.11 (CDS) was used for all data collection and processing. The power function value (PFV) was set in the program method conditions and stored in the audit trail. The response curves were generated at a minimum of 6 different PFVs, with most being run at 7. This resulted in the generation and processing of ~100 response curves.

Results

Linearity

All 14 data sets showed an increase in linear dynamic range. The PFV that returned the greatest correlation coefficient was then listed as the optimal value. Figure 2 shows the response curves for bovine serum albumin (BSA) sample analyzed using a monolith column chemistry on the CAD without a power transformation (PFV=1.0) and with a PFV of 1.7. The data without the PFV=1.0 was processed with a 2nd order polynomial fit (quadratic in CDS software), while the PFV=1.7 was processed with a standard linear regression. Neither used the origin as a data point. The results showed an improvement in correlation when using the power transformation over the 2.5 orders of magnitude analyzed.

FIGURE 2: Nine point response curves for bovine serum albumin.



The optimal PFV value and the resulting correlation coefficient for all samples analyzed over the full range are presented in Table 1A. The average PFV for these samples was 1.55. The values in green represented data points that had acceptable back calculations from the linear equation of the response curves. While the R^2 values for all 14 curves were greater than 0.995, many of the lower level back calculations were not within the 5% range and are colored according to the percent deviation. The highest PFV used was 1.7, which was not high enough for the pyridoxine sample resulting in poor back calculations. Overall, the samples showed a similar trend: good correlation for high mass concentrations but declining when decreasing by two or three orders of magnitude. In Table 1B, the same data are analyzed but excluding the highest two levels. Over this reduced mass range the average of the optimal PFVs was 1.38. The majority of back calculations now fell within the acceptable range and the correlation coefficients improved.

The linearity and back calculation results demonstrate some of the advantages of the power transformation, as well as some guidelines for use. The use of a large PFV can give the appearance of linearity over a wide dynamic range but it is always important to understand how each point is weighted over the entire range. Note, with any curve over two orders of magnitude, the points at the top of the curve will have a greater impact on the correlation coefficient than those at lower levels. This can result in high R² values but poor back calculations. If this function is intended to be used to measure low levels of analytes based on area percent, the quality of the curve must be evaluated. The major improvement noticed in this work was the ability to obtain linear response over a larger targeted range than previously available. If a curve still does not correlate after a power transformation has been used (i.e., the PFV value was either too high or too low), a second order polynomial fit can be used (4 data points or more required). In most cases, this will provide an improved fit over the use of the 2nd order polynomial alone (data not shown), with the added benefit of additional chromatographic improvements (see below).

Table 1: Results obtained at the optimal power function value over (A) a large range of 2 orders of magnitude or greater. (B) smaller ~ 1.5 orders of magnitude range. The color key to the correlation coefficient (R²) and back-calculated percent recovery is shown below.

1A	Prandione	Hydrocortisone	Ketoprofen	Warfarin	Progesterone	Pyridone	Procainamide Cls	Dibufenac	BSA	Dibufenac Cls	Euric Acid	Stearic Acid	Painic Acid	Myric Acid
Power Function	1.6	1.5	1.5	1.5	1.4	1.7	1.7	1.6	1.7	1.7	1.4	1.6	1.6	1.3
R2 Value	0.9992	0.9996	0.9990	0.9997	0.9996	0.9952	0.9999	0.9991	0.9993	0.9999	0.9996	0.9996	0.9992	0.9993
Std 1	99.9%	100.6%	100.9%	100.5%	99.4%	97.3%	100.0%	99.4%	98.9%	99.8%	100.1%	99.9%	99.5%	100.2%
Std 2	101.7%	98.1%	97.8%	98.9%	102.5%	107.8%	100.5%	103.9%	103.9%	99.8%	100.8%	101.5%	103.3%	98.7%
Std 3	97.2%	98.3%	95.0%	97.9%	101.2%	113.5%	99.7%	97.2%	102.5%	101.9%	96.8%	96.2%	98.8%	99.1%
Std 4	87.4%	95.5%	93.0%	93.8%	94.3%	#N/A	96.7%	91.8%	98.1%	99.2%	93.4%	95.5%	90.7%	108.7%
Std 5	91.2%	99.2%	98.0%	98.1%	88.2%	90.3%	93.8%	85.8%	94.3%	93.8%	93.7%	90.7%	84.4%	122.1%
Std 6	119.2%	113.3%	127.6%	114.5%	98.1%	65.2%	94.4%	96.9%	84.5%	88.1%	101.4%	95.7%	93.1%	112.1%
Std 7	192.0%	161.9%	200.1%	165.5%	134.5%	4.2%	115.2%	131.4%	77.4%	93.6%	137.3%	131.5%	133.0%	43.0%
Std 8	#N/A	#N/A	#N/A	#N/A	#N/A	-79.6%	174.3%	214.5%	88.5%	119.6%	223.0%	225.3%	234.6%	-59.3%
Std 9	#N/A	#N/A	#N/A	#N/A	#N/A	#N/A	#N/A	#N/A	#N/A	191.5%	#N/A	#N/A	#N/A	#N/A

1B	Prandione	Hydrocortisone	Ketoprofen	Warfarin	Progesterone	Pyridone	Procainamide	Dibufenac	BSA	Dibufenac Cls	Euric Acid	Stearic Acid	Painic Acid	Myric Acid
Power Function	1.2	1.3	1.2	1.2	1.2	1.7	1.6	1.3	1.7	1.6	1.2	1.4	1.3	1.5
R2 Value	0.9999	0.9999	0.9999	0.9997	0.9997	1.0000	0.9999	0.9999	0.9998	0.9999	0.9997	0.9994	0.9998	0.9988
Std 1	#N/A	#N/A	#N/A	#N/A	#N/A	#N/A	#N/A	#N/A	#N/A	#N/A	#N/A	#N/A	#N/A	#N/A
Std 2	#N/A	#N/A	#N/A	#N/A	#N/A	#N/A	#N/A	#N/A	#N/A	#N/A	#N/A	#N/A	#N/A	#N/A
Std 3	99.8%	100.0%	100.1%	99.5%	100.1%	100.0%	100.2%	99.8%	100.4%	100.1%	99.5%	99.4%	99.8%	99.7%
Std 4	100.8%	100.0%	100.3%	100.3%	100.3%	#N/A	99.6%	101.3%	98.5%	99.5%	102.3%	103.0%	101.3%	99.7%
Std 5	101.4%	97.5%	97.2%	100.9%	96.5%	98.9%	98.6%	98.9%	99.3%	100.4%	100.1%	98.5%	97.7%	106.8%
Std 6	97.0%	98.6%	99.7%	96.5%	97.8%	102.1%	97.4%	98.1%	99.0%	97.4%	95.3%	94.8%	97.0%	104.8%
Std 7	97.6%	107.8%	108.2%	93.9%	113.9%	99.6%	110.9%	101.8%	110.5%	105.6%	98.7%	101.6%	107.3%	70.5%
Std 8	#N/A	#N/A	#N/A	#N/A	#N/A	#N/A	#N/A	#N/A	#N/A	#N/A	#N/A	#N/A	#N/A	#N/A
Std 9	#N/A	#N/A	#N/A	#N/A	#N/A	#N/A	#N/A	#N/A	#N/A	#N/A	#N/A	#N/A	#N/A	#N/A

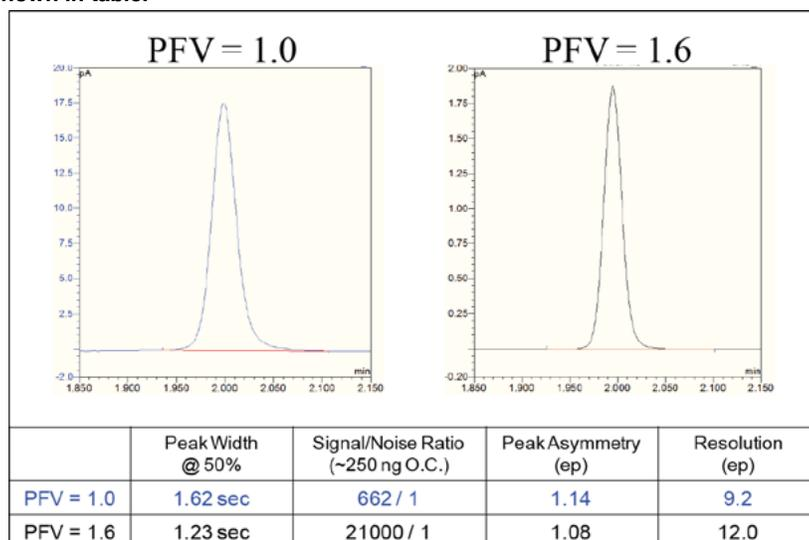
< 80%	80% - 90%	90% - 95%	95% - 105%	105% - 110%	110% - 120%	>120%
-------	-----------	-----------	------------	-------------	-------------	-------

Signal-to-Noise and Limits Testing

The impact of applying a power transformation to a number greater than 1 is to increase its value (e.g., with a PFV=1.3; 2 becomes 2.5, 20 becomes 49.1, and 200 becomes 280.2). However, the impact of applying that same power function to a value less than 1 is to decrease its value (e.g., 0.2 becomes 0.123). This mathematical principle along with the need to apply a scaling factor in order to keep all the data within the range of the detector, results in a dramatic decrease in the observed noise. While this has a greater impact on larger peaks, smaller peaks can also benefit. The degree to which the S/N was affected for the entire analytical range was analyzed.

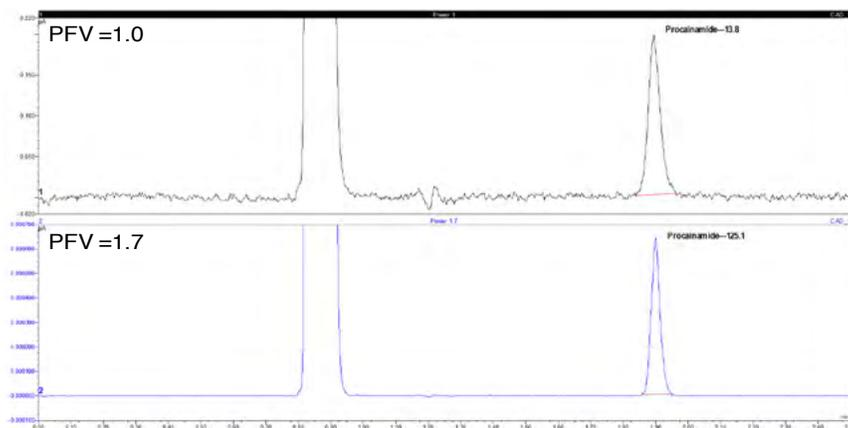
The signal-to-noise and peak width at both half height and base were measured and evaluated for every sample point at all the PFVs used above. This resulted in more than 750 S/N data points. For all 14 samples and across all mass ranges, the S/N was found to increase as the PFV increased. In general it was observed that the S/N was 1-2 times greater at the low end and 50-1000 times greater for high mass values (with a PFV of 1.5). However, the effect on S/N can vary depending on the baseline noise and gradient effects. For example, Figure 4 shows data for stearic acid analyzed as part of a mix of fatty acid standards using a water/acetonitrile to acetonitrile/IPA gradient. The S/N for the 250 ng on-column injection improved by a factor of 31 as shown.

FIGURE 4: Analysis of stearic acid with PFVs of 1.0 and 1.6 with peak attributes shown in table.



It is important to ensure that no value is lost by application of the power function. Peaks that are near or below the limit of detection (LOD) can be harder to discern after a power transformation. For this reason, trace analysis of unknown samples should initially be analyzed with a PFV=1.0 in order to ensure accuracy of the data. This however does not mean that the power transformation has no value at low levels. Figure 5 shows the analysis of procainamide at a PFV of 1.0 and 1.7. The S/N for the none-power transformed data of 13ng on column procainamide is just above the limit of quantification (LOQ) at 18.8/1 but with the PFV=1.7, the S/N was 126.1/1. This 13 ng on-column amount represents a 0.1% w/w impurity level for a 13 μ g API injection. Setting system suitability requirements for S/N at 20/1 (2x LOQ) for a point below the release criteria is needed for a robust method. In this example the 0.05 w/w% level would now meet the system suitability requirements and result in a more robust method.

FIGURE 5: Analysis of procainamide with power function values of 1.0 and 1.7 illustrating signal-to-noise improvements at low levels.

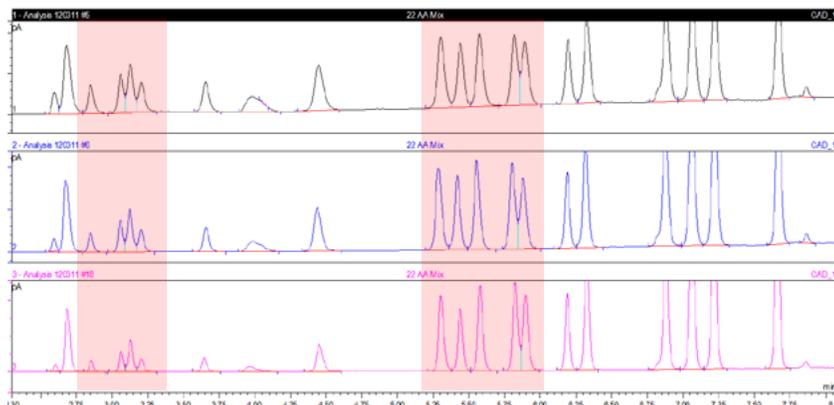


Peak Width and Resolution

The effect of the power transformation on other chromatographic results was also observed during the analysis of the linearity samples. The effect on peak shape also correlated as expected: greater impact on higher levels with the peaks becoming taller and narrower. This can be seen for the stearic acid peak (Figure 4), which showed a decreased peak width at half height of 24%. This narrowing has two effects on data quality. First, is an improved peak symmetry and second the narrowing of peaks throughout the chromatogram results in an improvement in the calculated resolution. The mathematical impact of power transformations on peak resolution is discussed in detail in work from Dasgupta.² This is experimentally shown again with the data from the fatty acid mix (Figure 4) but examined in a more complex-amino acid sample as well.

Figure 6 illustrates the analysis of 21 underivatized amino acids using the Corona ultra RS detector at PFVs of 1.0, 1.5 and 2.0. The two areas highlighted in red accounted for the regions with poorest peak resolution. Peak resolution was calculated by the software using the EP method with calculations based on next integrated peak. The resolution increased as the selected PFV increased with an average increase of 20% at PFV=1.5 and 39% at PFV=2.0. This improvement in the observed data for linearity, signal-to-noise, and resolution can be implemented without negatively impacting data quality or accuracy for these samples.

FIGURE 6: Analysis of 21 underivatized amino acids using reversed phase chemistry from 0.4% heptafluorobutyric acid to 50% acetonitrile at multiple PFVs. The peak resolution was calculated for all peaks highlighted in red and the results are shown in table below.



Resolution of Selected Peaks						
	2.85 min	3.06 min	5.30 min	5.42 min	5.57 min	5.81 min
PFV = 1.0	2.99	0.88	1.62	1.61	2.77	0.78
PFV = 1.5	3.64	1.07	1.81	2	3.39	0.96
PFV = 2.0	4.15	1.28	2.17	2.28	3.79	1.11

Conclusion

- Implementation of a power transformation to raw data with charged aerosol detection results in improvements in observed linear range, signal-to-noise, and peak resolution, along with reduction in peak width and asymmetry.
- The power function value should be chosen based on the linear range needed and experimental results as one preset value is not suited for all analyses.
- The improvement in signal-to-noise at levels above the limit of quantification provides a powerful tool for improving reproducibility and quality when doing limits testing with the Corona ultra RS charged aerosol detector.
- The data presented here illustrated that use of power transformation in the Corona ultra RS is auditable and can improve data quality.

References

1. Górecki, T.; Lynen, F.; Szucs, R.; Sandra, P. Universal Response in Liquid Chromatography Using Charged Aerosol Detection. *Anal. Chem.* **2006**, *78*, 3186–3192.
2. Dasgupta PK; Chen Y; Serrano CA; Guiochon G; Liu H; Fairchild JN; Shalliker RA. Black Box Linearization for Greater Linear Dynamic Range: The Effect of Power Transforms on the Representation of Data. *Anal Chem.* **2010** Dec 15;82(24):10143-50.
3. Shalliker RA; Stevenson PG; Shock D; Mnatsakanyan M; Dasgupta PK; Guiochon G. Application of Power Functions to Chromatographic Data for the Enhancement of Signal to Noise Ratios and Separation Resolution. *J. Chrom A*, **2010**, 1217, 5693-5699.

www.thermofisher.com/dionex

©2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.

U.S./Canada	(847) 295 7500	Denmark	(45) 36 36 90 90	Sweden	(46) 8 473 3380	India	(91) 22 2764 2735
Brazil	(55) 11 3731 5140	France	(33) 1 39 30 01 10	Switzerland	(41) 62 205 9966	Japan	(81) 6 6885 1213
Austria	(43) 1 616 51 25	Germany	(49) 6126 991 0	United Kingdom	(44) 1276 691722	Korea	(82) 2 2653 2580
Benelux	(31) 20 683 9768	Ireland	(353) 1 644 0064	Australia	(61) 2 9420 5233	Singapore	(65) 6289 1190
	(32) 3 353 42 94	Italy	(39) 02 51 62 1267	China	(852) 2428 3282	Taiwan	(886) 2 8751 6655

Thermo
S C I E N T I F I C

Part of Thermo Fisher Scientific

PN70003_E 07/16S