

Industrial

Automated, multi-parameter gasoline characterization using GC-VUV and ASTM D8071

Authors

Daniela Cavagnino¹, Oliver Reyes²,
Amit Gujar², Peter Boler³, Sean Jameson³

¹Thermo Fisher Scientific, Milan, Italy

²Thermo Fisher Scientific, Texas, USA

³VUV Analytics, Cedar Park, Texas, USA

Keywords

Gas chromatography, VUV spectroscopy, PIONA analysis, hydrocarbons, ASTM D8071, fuel composition, paraffins, olefins, aromatics, ethanol, benzene, GC-VUV, D1319, D5599, D5769, D3606, D6550

Goal

This application note details an innovative method for fully automated PIONA class analysis and hydrocarbon profiling of gasoline using gas chromatography with vacuum ultraviolet detection (GC-VUV) under ASTM D8071.¹ The method delivers comprehensive compositional data—including paraffins, isoparaffins, olefins, naphthenes, aromatics, ethanol, and benzene—in a single 34-minute run. With no sample prep, calibration standards, or multiple instruments required, this platform significantly reduces cost-per-sample and complexity, while improving precision and compliance with global fuel regulations.

Introduction

Gasoline is a complex hydrocarbon blend whose composition directly impacts combustion behavior, emissions, and regulatory compliance. Accurately quantifying components such as aromatics, olefins, ethanol, and benzene is critical, but traditional ASTM methods require multiple techniques (FIA, GC-FID, GC-MS, etc.) to measure all relevant parameters. Table 1 summarizes several test methods and their respective scopes.

ASTM D8071 introduces a consolidated method using GC-VUV, offering improved precision, spectral resolution, and reduced operational burden. This approach is now recognized as an EPA- and CARB-approved alternative to several legacy test methods, including ASTM D1319, D5599, D5769, and D3606.

Table 1. Several gasoline test methods and their parameters

| ASTM method | D6550 | D4815 | D5599 | D1319 | D3606 | D5769 | D5580 | D6729 / D6730 | D6839 | D8071 |
|--------------------|-------|----------|---------|-------|--------|-------|----------|---------------|---------------|--------|
| Technique | SFC | MDGC-FID | GC-OFID | FIA | GC-TCD | GC-MS | MDGC-FID | GC-FID | Reformulyzer™ | GC-VUV |
| Aromatics | | | | ✓ | | ✓ | ✓ | ✓ | ✓ | ✓ |
| Benzene | | | | | ✓ | ✓ | ✓ | ✓ | ✓ | ✓ |
| Olefins | ✓ | | | ✓ | | | | ✓ | ✓ | ✓ |
| Ethanol | | ✓ | ✓ | | | | | ✓ | ✓ | ✓ |
| Ethyl Benzene | | | | | | ✓ | ✓ | ✓ | | ✓ |
| Isoparaffin | | | | | | | | ✓ | ✓ | ✓ |
| Methanol | | ✓ | ✓ | | | | | ✓ | | ✓ |
| Methyl Naphthalene | | | | | | | | ✓ | | ✓ |
| Naphthalene | | | | | | | | ✓ | | ✓ |
| Naphthene | | | | | | | | ✓ | ✓ | ✓ |
| Paraffin | | | | | | | | ✓ | ✓ | ✓ |
| Toluene | | | | | ✓ | ✓ | ✓ | ✓ | | ✓ |

This application note describes an approach to achieving PIONA class and multi-parameter analysis with a single measurement using the VUV Analyzer Platform for Fuels running ASTM D8071.²

Experimental

Gasoline samples were analyzed on the VUV Analyzer Platform for Fuels with no sample preparation required. The VUV Analyzer consists of the Thermo Scientific™ TRACE™ 1610 gas chromatograph coupled to the VGA-100™ spectrometer detector (VUV Analytics, Inc.).

Data are acquired and processed with the VUVision™ Software and VUV Analyze™ Software (VUV Analytics, Inc), respectively, specifically configured to run ASTM D8071.

Workflow overview

The VUV Analyzer Platform for Fuels performs gasoline PIONA class analysis using a streamlined five-step workflow:

- 1. System validation** – Verifies instrument performance using standard hydrocarbons.
- 2. Sample prep** – No preparation required.
- 3. Data acquisition** – Performed using VUVision Software.
- 4. Spectral matching** – Automated deconvolution with VUV Analyze Software.
- 5. Quantitation** – Volume % and mass % calculated with high accuracy using built-in response factors and density data.

Results and discussion

Three-dimensional spectral analysis

Traditional chromatography identifies and quantifies compounds using peak retention time and peak tables. Because of this, it is important that peaks of interest are sufficiently baseline resolved. As gasoline is a complex mixture of hydrocarbons, achieving sufficient baseline resolution can be difficult as numerous compounds tend to coelute. Using a longer column can provide better separation, but it will also extend run times. Alternatively, other techniques sometimes require the use of complex valves, multiple columns, and compound and class-specific traps.

Unlike conventional GC, GC-VUV collects data across three axes: retention time, absorbance, and wavelength. Every compound has a unique spectral fingerprint, and members of the same hydrocarbon class (e.g., aromatics or olefins) exhibit similar absorbance patterns. This spectral distinction allows for accurate class separation even in the presence of coelution. As a result, analysis of gasoline can be accomplished quickly using a single 30-meter column solution.

Reformulated gasoline samples containing over 300 compounds were analyzed using a single-column setup in just 34 minutes, as shown in Figure 1.

Class-specific spectral filters are applied to overlay the chromatogram and distinguish paraffins, isoparaffins, olefins, naphthenes, and aromatics visually and analytically. Class-based spectra can be combined to provide accurate class-based analysis that is required for PIONA.

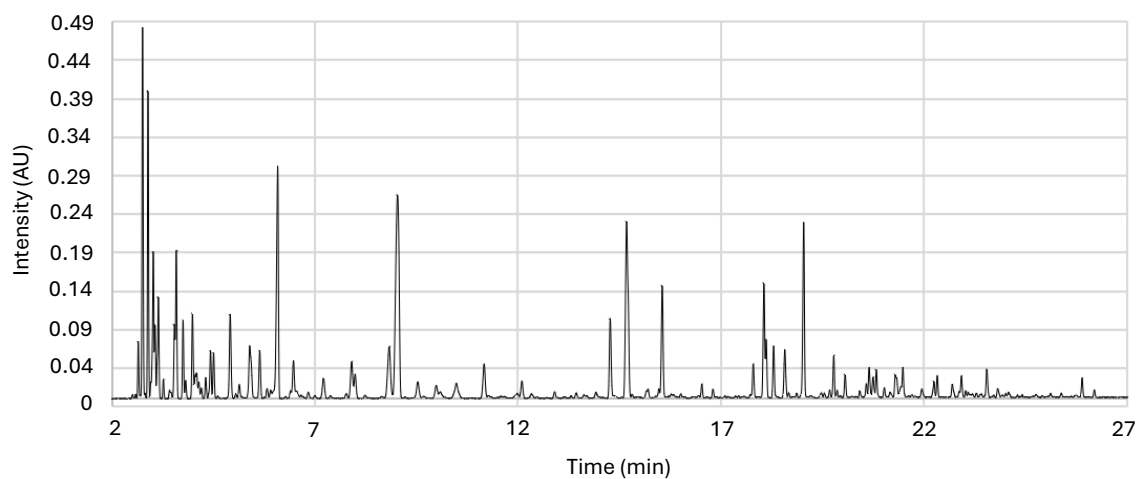


Figure 1. Reformulated gasoline sample with no spectral filters applied

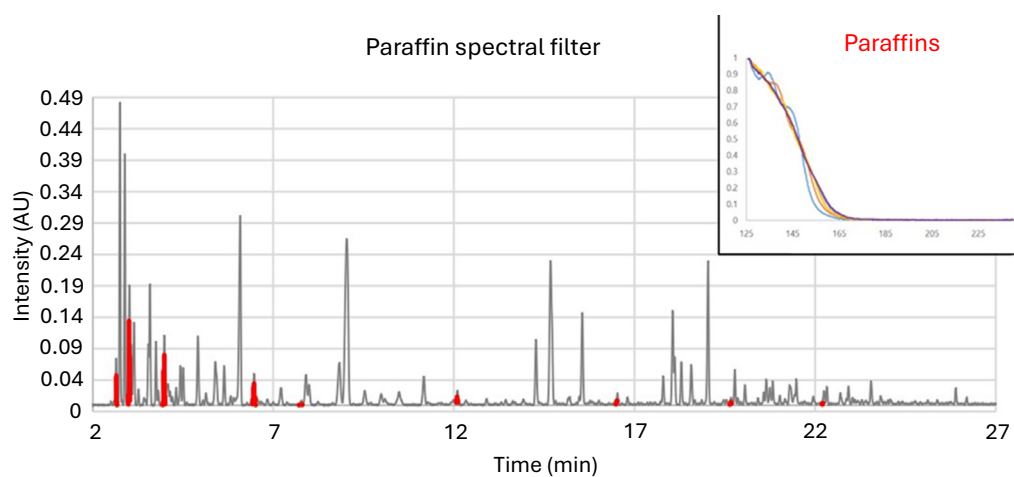


Figure 2. Reformulated gasoline chromatogram with paraffin spectral filter applied. The inset shows the VUV absorbance spectra of several common paraffins.

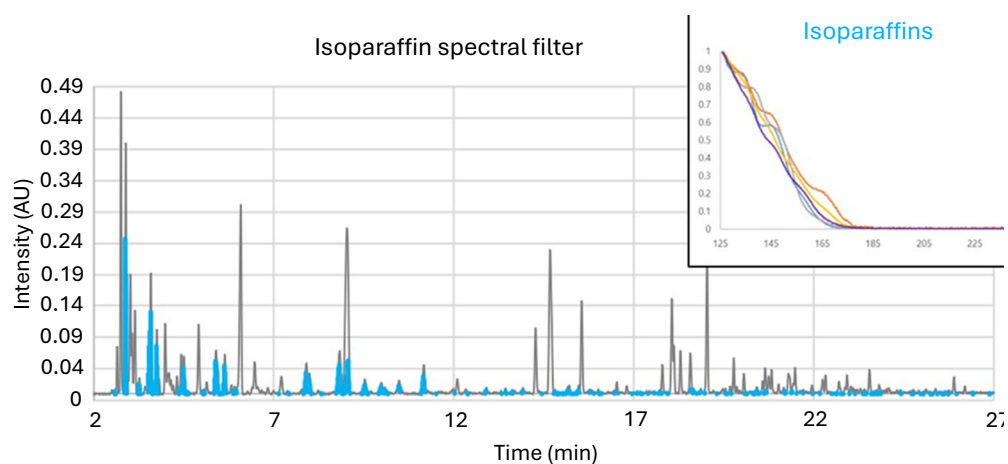


Figure 3. Reformulated gasoline chromatogram with isoparaffin spectral filter applied. The inset shows the VUV absorbance spectra of several common isoparaffins.

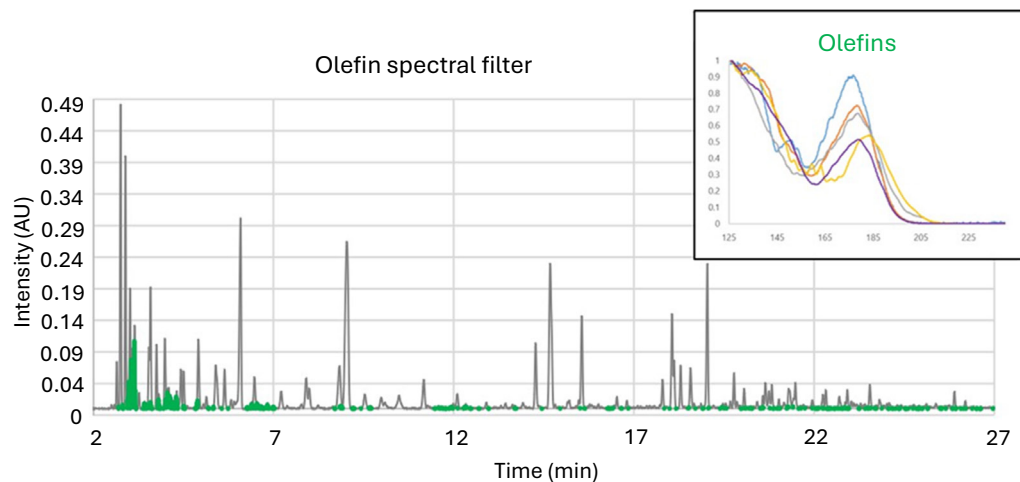


Figure 4. Reformulated gasoline chromatogram with olefin spectral filter applied. The inset shows the VUV absorbance spectra of several common olefins.

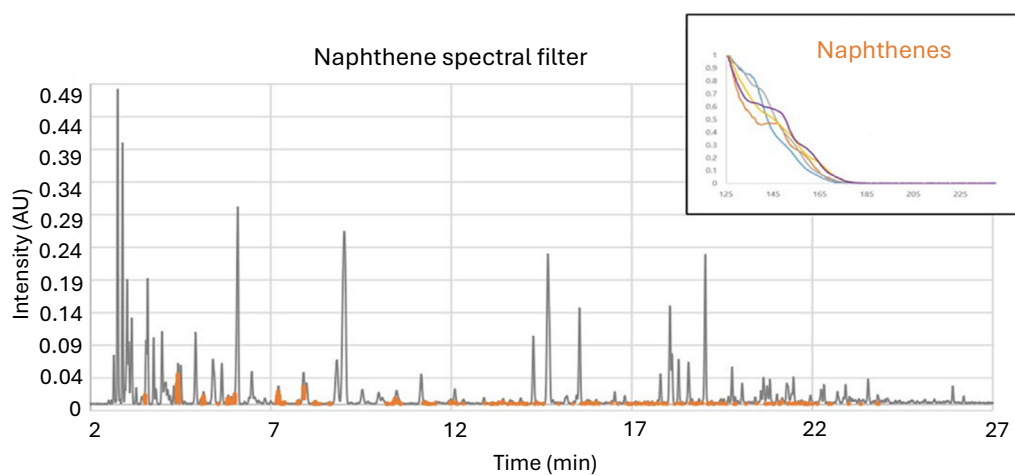


Figure 5. Reformulated gasoline chromatogram with naphthene spectral filter applied. The inset shows the VUV absorbance spectra of several common naphthenes.

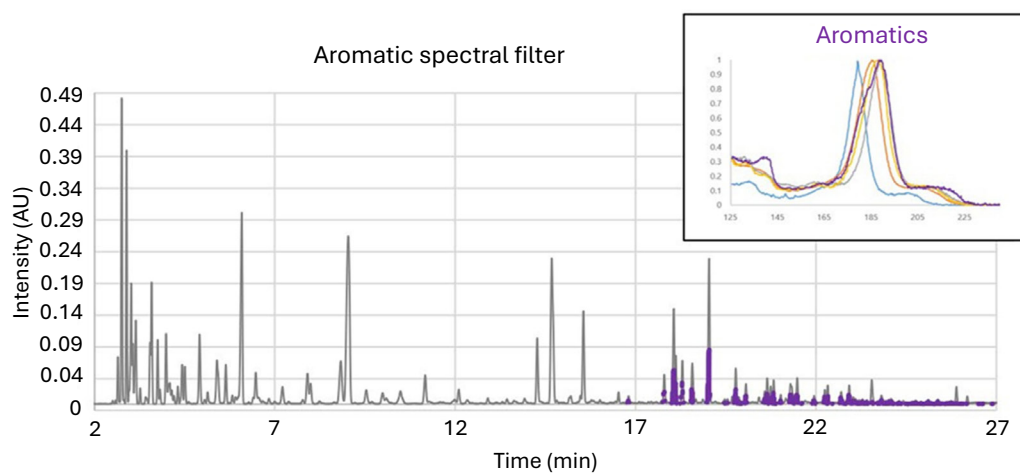


Figure 6. Reformulated gasoline chromatogram with aromatic spectral filter applied. The inset shows the VUV absorbance spectra of several common aromatics.

Figures 2–6 display the spectral filters associated with each of the PIONA classes along with an overlay of the individual spectra that are used in that filter. As one can see, spectra of a given class share similar shapes. Spectral filters give a good representation of where and when compounds of a given class absorb in the GC-VUV chromatogram.

Several classes of compounds coelute in the GC-VUV chromatogram. However, with the VUV Analyzer Platform, the GC-VUV chromatogram is divided into regularly spaced time intervals during analysis. Each spectrum can then be automatically compared and matched against a compound library and analyzed to determine the contribution of each compound. This automated approach is called Time Interval Deconvolution™, and it allows for accurate class-based analysis. When a coelution occurs, VUV Analyze Software uses the unique spectral shapes of each class and compound to determine the best multi-analyte fit.

After VUV Analyze Software completes the carbon number and class categorization of the components within the sample, an automated calculation determines the mass percent and volume

percent makeup. The result is a table based on compound class (Table 2).

Individual compound speciation

In addition to class-based reporting, the VUV Analyzer Platform running ASTM D8071 provides detailed insight into key compounds of interest. For fuels certification, those compounds include methanol, ethanol, benzene, iso-octane, toluene, ethylbenzene, naphthalene, methylnaphthalene, and xylenes (Table 3), with the option to add other compounds for non-regulated purposes. Individual speciated components can be identified, even if they coelute, using the Time Interval Deconvolution described above.

Key outcomes

- Accurate quantification of PIONA classes with coelution resolved via spectral deconvolution
- Ethanol, benzene, toluene, xylenes, and other regulatory targets detected and quantified
- Results presented as both volume % and mass % across carbon number
- Easy-to-read tables and overlays simplify data interpretation

Table 2. ASTM D8071 results are presented in an easy-to-read table showing carbon number and class breakdown. Note that ethanol is not included in the PIONA table. As a result, totals equal only 88.9.

| Mass | | | | | | |
|--------|--------|---------|--------|--------|---------|---------|
| C. no. | P | I | O | N | A | Total |
| C1 | | | | | | |
| C2 | | | | | | |
| C3 | | | | | | |
| C4 | 1.1144 | 0.1064 | 0.0450 | | | 1.2657 |
| C5 | 2.9113 | 6.0742 | 3.7817 | 0.2446 | | 13.011 |
| C6 | 2.0138 | 6.9951 | 2.2886 | 1.6600 | 0.8034 | 13.7610 |
| C7 | 1.2529 | 5.5536 | 0.8038 | 2.0491 | 3.8278 | 13.4871 |
| C8 | 0.5290 | 17.2473 | 0.7076 | 2.0670 | 5.7730 | 26.3239 |
| C9 | 0.3473 | 3.2496 | 0.1982 | 0.9024 | 5.9828 | 10.6802 |
| C10 | 0.2080 | 0.8515 | 0.6075 | 0.9567 | 3.1507 | 5.7743 |
| C11 | 0.1045 | 0.9554 | 0.3179 | 0.5486 | 1.3512 | 3.2776 |
| C12 | | 0.1992 | 0.1638 | 0.0571 | 0.3897 | 0.8098 |
| C13 | | 0.1849 | 0.0491 | | 0.2292 | 0.4633 |
| C14 | | 0.0445 | | | | 0.044 |
| C15 | | 0.0028 | | | | 0.0028 |
| C16 | | | | | | |
| C17 | | | | | | |
| C18 | | | | | | |
| C19 | | | | | | |
| Total | 8.4811 | 41.4645 | 8.9633 | 8.4855 | 21.5077 | |

| Volume % | | | | | | |
|----------|--------|---------|--------|--------|---------|---------|
| C. no. | P | I | O | N | A | Total |
| C1 | | | | | | |
| C2 | | | | | | |
| C3 | | | | | | |
| C4 | 1.4101 | 0.1398 | 0.0538 | | | 1.6037 |
| C5 | 3.4051 | 7.1802 | 4.1864 | 0.2404 | | 15.0121 |
| C6 | 2.2368 | 7.7921 | 2.3616 | 1.6090 | 0.6695 | 14.6690 |
| C7 | 1.3422 | 5.9535 | 0.8039 | 1.9908 | 3.2336 | 13.3239 |
| C8 | 0.5515 | 17.9972 | 0.7041 | 1.9749 | 4.8632 | 26.0909 |
| C9 | 0.3545 | 3.3225 | 0.1979 | 0.8451 | 5.0351 | 9.7550 |
| C10 | 0.2087 | 0.8459 | 0.5969 | 0.8663 | 2.6164 | 5.1341 |
| C11 | 0.1034 | 0.9195 | 0.3101 | 0.4966 | 1.1072 | 2.9368 |
| C12 | | 0.1892 | 0.1581 | 0.0511 | 0.3162 | 0.714 |
| C13 | | 0.1741 | 0.0469 | | 0.1925 | 0.4135 |
| C14 | | 0.0416 | | | | 0.0416 |
| C15 | | 0.0026 | | | | 0.0026 |
| C16 | | | | | | |
| C17 | | | | | | |
| C18 | | | | | | |
| C19 | | | | | | |
| Total | 9.6122 | 44.5583 | 9.4195 | 8.0742 | 18.0337 | |

Table 3. Individual speciated compounds identified in reformulated gasoline sample

| Report item | Category | Retention time (min) | Mass% | Volume % |
|--------------------|-------------|----------------------|---------|----------|
| Methanol | Alcohol | - | - | - |
| Ethanol | Alcohol | 2.7468 | 11.0980 | 10.3021 |
| Benzene | Aromatic | 4.9224 | 0.8034 | 0.6695 |
| Iso-octane | Isoparaffin | 6.0768 | 7.2400 | 7.6640 |
| Toluene | Aromatic | 9.0894 | 3.8278 | 3.2336 |
| Ethylbenzene | Aromatic | 14.2686 | 1.0800 | 0.9123 |
| Naphthalene | Aromatic | 23.5260 | 0.2053 | 0.1467 |
| Methylnaphthalenes | Aromatic | - | 0.2372 | 0.1725 |
| Xylenes | Aromatic | - | 4.6930 | 3.9508 |

Table 4. The applicable test ranges for ASTM D6708 correlation equations

| Correlated method | Units | ASTM D8071 reported results | | |
|------------------------|----------|-----------------------------|--------|--------------------------------|
| | | Min | Max | Correlation equation |
| ASTM D1319 (aromatics) | % Volume | 14.743 | 58.124 | (D8071 vol%) - 0.8313 |
| ASTM D1319 (olefins) | % Volume | 0.0190 | 17.412 | (D8071 vol%) x 0.827 + 0.5318 |
| ASTM D3606 (benzene) | % Volume | 0.12 | 0.946 | (D8071 vol%) x 0.968 - 0.016 |
| ASTM D5599 (ethanol) | Mass % | 0.396 | 15.991 | (D8071 mass%) - 0.0878 |
| ASTM D5769 (aromatics) | % Volume | 14.743 | 36.068 | (D8071 vol%) - 1.097 |
| ASTM D5769 (benzene) | % Volume | 0.09 | 1.09 | (D8071 vol%) x 0.9446 - 0.0014 |
| ASTM D6550 (olefins) | Mass % | 0.24 | 16.71 | (D8071 mass%) + 0.2979 |

The applicable ASTM D8071 test result range for ASTM D6708 correlation equations, as shown in Table 4, provide equivalence to traditional referee methods. Regulatory agencies, including the EPA, CGSB, and CARB, have accepted this method as an alternative for multiple compound classes.

Cost advantages

Cost-per-analysis using GC-VUV running ASTM D8071 is significantly less on a per-sample basis than the alternative approaches. This is due in large part to lower labor and consumables costs resulting from the elimination of sample preparation and complex apparatus setup, automation provided by the VUV Analyze Software, minimal ongoing consumable and maintenance costs, and the consolidation of multiple techniques into a single, easy-to-use method.

To better compare cost-per-analysis, ASTM D8071 was compared to ASTM D1319, ASTM D3606, ASTM D5599, and ASTM D5769, each of which ASTM D8071 has ASTM D6708 equivalency with. To ensure consistency, the parameters used in the calculations included: capital cost of the analytical hardware depreciated over five (5) years, a consistent utilization rate of 80% across all techniques, expected annual consumable and maintenance costs across all techniques, and the cost of labor and labor time spent per day interfacing with a given technique.

Figure 7 outlines the results of that analysis and shows a substantial difference in cost-per-analysis between ASTM D8071 and the alternative methodologies required to acquire the same data set. While having similar capital costs, the consumables and labor costs when using ASTM D8071 are over 12 times less expensive to run on a per-sample basis compared to the alternative methodologies.

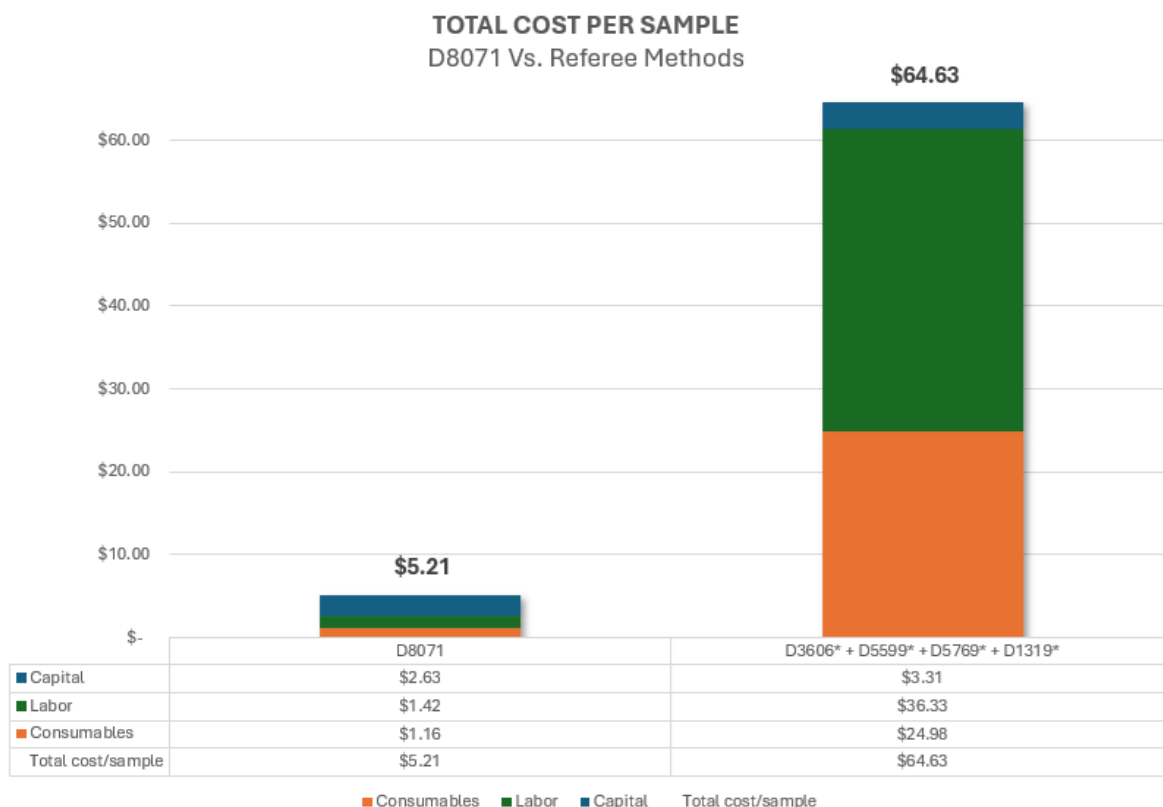


Figure 7. Cost-per-sample analysis. Note that all values are in USD.

Conclusion

GC-VUV with ASTM D8071 enables complete hydrocarbon group and compound-level analysis of gasoline in a single method. Its ability to automate spectral deconvolution and report mass and volume percentages simplifies operations while meeting strict compliance standards. This solution not only enhances analytical accuracy and repeatability but also drastically reduces operating costs and lab complexity.

- Short analysis time with full gasoline class and compound profiling in 34 minutes
- Faster throughput with full automation from injection to reporting
- 12× lower cost-per-analysis, due to reduced labor, no calibration standards, and consolidated workflow

References

1. ASTM International - D8071 Standard Test Method
2. VUV Analytics – ASTM D8071 for Gasoline Analysis

Learn more at thermofisher.com/gc-vuv