

Thermo Scientific™ Prima PRO Process Mass Spectrometer

Fast On-line monitoring of flare gases

Author: Graham Lewis, Thermo Fisher Scientific, Winsford, Cheshire, United Kingdom

Keywords

- Lower Heating Value (LHV)
- Net Heating Value
- Higher Heating Value (HHV)
- Gross Heating Value
- Environmental Protection Agency (EPA)
- EPA 40 CFR Parts 60 and 63
- EPA Refinery Sector Rule (RSR) 40 CFR 63
- Magnetic sector
- Rapid multistream sampler
- Total sulfur

Introduction

Increasingly, national and local government bodies are requiring processing companies to monitor the various emissions from their plant stacks and flares to reduce the amount of pollution entering the atmosphere. For example, the US Environmental Protection Agency (EPA) has published amendments to Refinery Sector Rule (RSR) 40 CFR Part 63 affecting flares; refineries must bring flares into compliance with new §63.670 'Requirements for Flare Control Devices' by January 30, 2019. The new requirements define five flare operating limits—combustion zone net heating value (NHV_{CZ}), dilution net heating value (NHV_{dil}), flare tip exit velocity, pilot flame presence, and visible emissions—and specifies an NHV_{CZ} minimum operating limit of 270 BTU/scf, based on a 15 minute block period¹. NHV_{CZ} can be calculated by measuring the net heating value of the vent gas (NHV_{VG}), making flare gas analysis a vital part of any compliance strategy. If the NHV_{CZ} approaches 270 BTU/scf, additional fuel gas such



as propane or natural gas have to be added. This may then require the addition of steam to the flare, to avoid the production of visible emissions. The analysis of flare gases presents a series of challenges to the processing industries. Firstly, the emissions are typically made up of complex mixtures of inorganic and organic species, and compositions and concentrations vary dramatically over time as process conditions change. And although many regulations require just the total heating value, total sulfur, or total hydrocarbon values to be recorded, measuring the concentrations of individual components helps identify the source of the emission, locating the problem to a specific part of the plant. Finally, the analysis speed is crucial as the heating value of the flare can quickly change. Analysis times measured in minutes increase the risk of failing to meet

emission standards.

Measurement of flare gas streams by process mass spectrometry

Process Mass Spectrometry is particularly suited to the measurement of flare gas streams because it offers accurate, fast, multicomponent analysis. Table 1 shows an example of a flare gas stream containing hydrogen, nitrogen and hydrocarbons up to C₆. Analysis of these eleven components will typically be performed in just 20 seconds, allowing one mass spectrometer to monitor

more than one flare, depending on the distances involved. Alternatively, one MS can be used to measure process streams and flare streams, even though the composition and concentrations may be very different.

Advantages of Prima PRO Process MS

At the heart of the Prima PRO is a magnetic sector analyzer (Figure 1) which offers unrivalled precision and accuracy compared with other mass spectrometers. Thermo Fisher Scientific manufactures both quadrupole and magnetic

Table 1 Example of flare gas stream composition

Component			Flare gas Typical composition %mol
Name	Formula	Molecular weight	
Hydrogen	H ₂	2	1 - 3
Nitrogen	N ₂	28	40 - 85
Ethylene	C ₂ H ₄	28	5 - 20
Ethane	C ₂ H ₆	30	5 - 20
1-Butene	C ₄ H ₈	56	0.5 - 1.5
i-Butane	i-C ₄ H ₁₀	58	0.1 - 0.5
n-Butane	i-C ₄ H ₁₀	58	0.1 - 0.3
1-Hexene	C ₆ H ₁₂	84	2.5 - 7.5
n-Hexane	n-C ₆ H ₁₄	86	0 - 1

sector mass spectrometers; over thirty years of industrial experience have shown the magnetic sector based analyzer offers the best performance for industrial online gas analysis.

Key advantages of magnetic sector analyzers include improved precision, accuracy, long intervals between calibrations and resistance to contamination. Typically, analytical precision is between 2 and 10 times better than a quadrupole analyzer, depending on the gases analyzed and complexity of the mixture.

Neutral gas atoms and molecules are first converted into positively charged ions in the Prima PRO ion source. This is an enclosed type for high sensitivity, minimum background interference and maximum contamination resistance. It is a high-energy (1000 eV) analyzer that offers extremely rugged performance in the presence of gases and vapors that have the potential for contaminating the internal vacuum components. Prima PRO has a proven track record of monitoring high percent level concentrations of organic compounds without experiencing drift or contamination.

Ions are then accelerated through a flight tube, where they are separated by their mass to charge ratios in a magnetic field of variable strength. Since the magnetic

sector mass spectrometer produces a focused ion beam at the detector, the peak shape obtained is 'flat-topped' and uniform response is observed over a finite mass width. As the height of the peak is directly proportional to the number of ions striking the detector it is also directly proportional to the concentration of the component being measured. As long as the measurement is taken anywhere on the peak's flat top, high precision analysis will be observed.

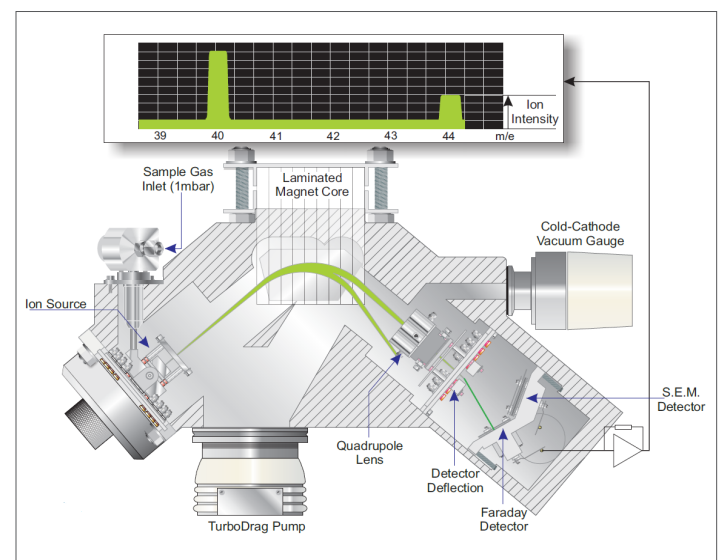


Figure 1 Prima PRO's magnetic sector analyzer

Prima PRO's ability to measure over a wide dynamic range is critically important if the varying composition levels in flare gas are to be measured accurately. Prima PRO has been independently evaluated by EfecTech UK, an independent specialist company providing accredited calibration and testing services to the energy and power

industries for gas quality, flow and total energy metering. It is accredited to internationally recognised ISO/IEC 17025:2005 standards; this specifies the general requirements for the competence to carry out tests and/or calibrations, including sampling.

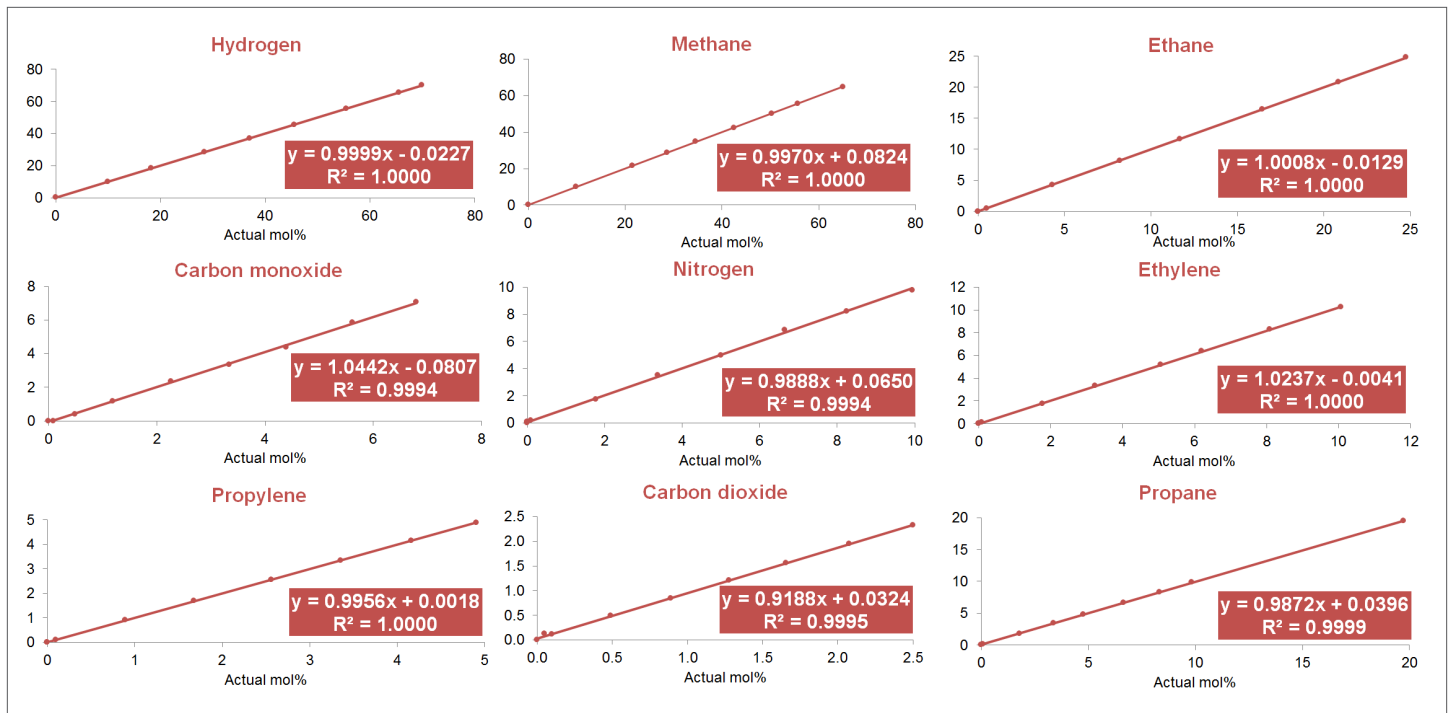


Figure 2 Prima PRO linearity data

Effectech linearity tests

Prima PRO was calibrated for sensitivity using just one gas mixture, and then tested with a set of nine reference gases covering a wide range of compositions. Figure 2 shows the linearity plots generated by EfecTech. They demonstrate significantly better linearity than that achieved by a thermal conductivity detector fitted to a gas chromatograph, and prove that Prima PRO is capable of generating accurate, reliable composition data from complex gas mixtures.

Table 2 Statistical report on Prima PRO 30 day run analyzing 13-component calibration cylinder without calibration

	Average	Standard Deviation	Relative Standard Deviation
Hydrogen	18.98	0.106	0.56%
Methane	15.65	0.068	0.44%
Acetylene	1.19	0.007	0.58%
Ethylene	33.77	0.072	0.21%
Ethane	10.09	0.054	0.53%
Methyl acetylene	0.39	0.008	2.12%
Propylene	5.73	0.014	0.24%
Carbon dioxide	0.44	0.017	3.72%
Propane	5.87	0.012	0.21%
Butadiene	0.99	0.006	0.59%
n-Butene	0.56	0.004	0.69%
i-Butene	0.50	0.012	2.50%
n-Butane	5.83	0.022	0.37%

Long-term stability

Prima PRO's long-term stability means the system can run for long periods without recalibration, significantly longer than quadrupole MS systems. Table 2 shows standard deviations achieved on a 30-day run without recalibration on a gas cylinder containing 13 different compounds, including 11 hydrocarbons.

Rapid multistream sampling

If the MS is to monitor multiple flare streams, or flare streams and process streams, then a fast, reliable means of switching between streams is required. Solenoid valve manifolds have too much dead volume and rotary valves suffer from poor reliability so we developed the unique Rapid Multistream Sampler (RMS). It offers an unmatched combination of sampling speed and reliability and allows sample selection from 1 of 32 or 1 of 64 streams. Stream

settling times are application dependent and completely user configurable. The RMS includes digital sample flow recording for every selected stream. This can be used to trigger an alarm if the sample flow drops—if a filter in the sample conditioning system becomes blocked, for example.

The RMS can be heated to 120 °C and the position of the stream selector is optically encoded for reliable, software controlled stream selection. Temperature and position control signals are communicated via Prima PRO's internal network.

The RMS has a proven track record of carrying out 10 million operations between maintenance and has a three year warranty as standard. No other multistream sampling device offers the same level of guaranteed reliability.

Measuring fuel properties with Prima PRO

The following fuel properties can routinely be derived by the Prima PRO mass spectrometer:

Lower Heating Value (LHV)	Also known as Net Heating Value or Lower Calorific Value
Higher Heating Value (HHV)	Also known as Gross Heating Value or Higher Calorific Value
Compressibility	
Actual Lower Heating Value	$\frac{\text{Ideal Lower Heating Value}}{\text{Compressibility}}$
Density	
Specific Gravity	
Lower Wobbe Index	$\frac{\text{Lower Heating Value}}{\sqrt{(\text{Specific Gravity})}}$
Higher Wobbe Index	$\frac{\text{Higher Heating Value}}{\sqrt{(\text{Specific Gravity})}}$
Air Requirement	
Combustion Air Requirement Index (CARI)	$\frac{\text{Air Requirement}}{\sqrt{(\text{Specific Gravity})}}$

Lower Heating Value is calculated using the value of 1212 BTU/scf for Hydrogen as recommended by EPA 40 CFR Parts 60 and 63, instead of the theoretical value of 274 BTU/scf. This provides a better indication of flare performance (e.g. meeting the minimum operating limit of 270 BTU/scf). Accurate analysis of hydrogen over a wide dynamic range is critical because of its high combustion potential; Prima PRO's magnetic sector analyzer is ideally suited to hydrogen analysis because it does not suffer from the 'zero blast' that makes analysis of light molecules problematic on many quadrupole analyzers.

HHV differs from LHV in that it takes into account the latent heat of vaporization of water in the combustion products.

The precision of these measurements with Prima PRO is normally better than 0.1% relative.

Figure 3 shows how the Ideal LHV of the flare gas stream is calculated by GasWorks® Derived Values, based on the individual components' Ideal LHVs.

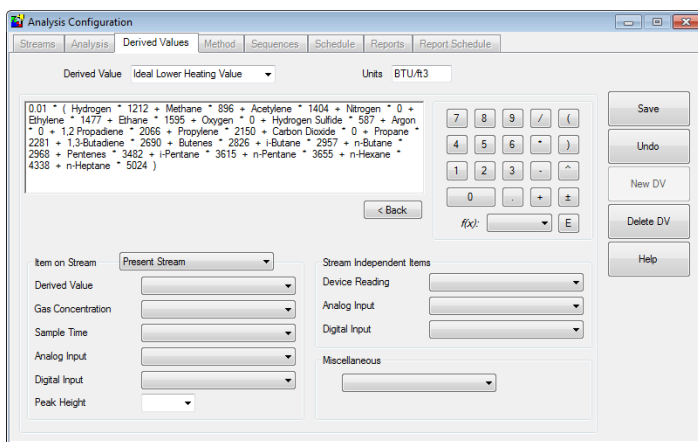


Figure 3 GasWorks® derived value for ideal lower heating value

Analysis times are typically 30 seconds or less, including settling time. Data are communicated to the plant host computer as they are measured, by one or more of a number of available methods, e.g. 4-20 mA or 0-10V analog outputs, Modbus, Profibus, or OPC.

Analytical Set-up

Thermo Scientific GasWorks software supports an unlimited number of analysis methods, enabling the analysis to be optimized on a per-stream basis. The most efficient peak measurements and the most appropriate speed versus precision settings can be selected for each gas stream, depending on process control requirements. Examples of different analysis methods are shown in Figures 4 and 5. Figure 4 shows an example of the analytical method for a flare gas stream, Figure 5 shows that for a process stream. Both methods, along with additional process methods, are being utilized on a single Prima PRO process MS measuring a combination of process streams and flare gas.

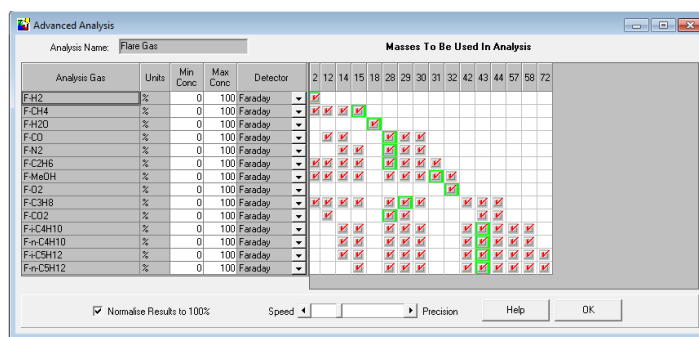


Figure 4 Example of flare stream method

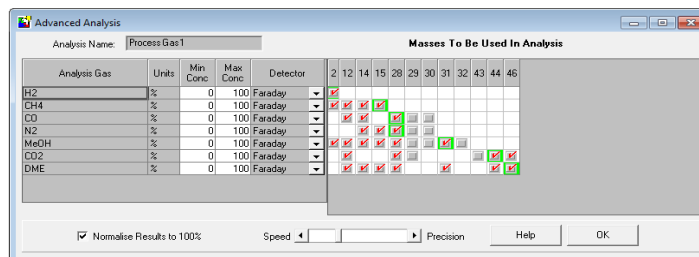


Figure 5 Example of process stream method

These two methods clearly show the amount of spectral overlap in the MS fragmentation patterns. It is extremely important to measure these fragmentation patterns from the actual components of interest—the use of surrogate compounds may simplify the calibration process but will inevitably lead to a reduction in accuracy.'

Test Data & Process Data

Table 3 shows Prima PRO's analytical performance on the two different stream types displayed in Figures 4 and 5. Stream analysis cycle time was 30 seconds. If multiple sample streams are measured the stream switching time and settling time is 10 seconds, so n streams can be analyzed with a cycle time of n x 40 seconds. Analytical performance was demonstrated during commissioning by analyzing two calibration gases periodically over 8 hours and observing the standard deviation values detailed in the table.

Table 3 Example of Prima PRO performance specification for flare gas and process gas streams

Component	Formula	Flare gas		Process gas	
		Concentration %mol	Prima PRO precision (standard deviation) %mol ≤	Concentration %mol	Prima PRO precision (standard deviation) %mol ≤
Hydrogen	H ₂	Balance	0.1	2	0.01
Methane	CH ₄	10	0.02	Balance	0.05
Carbon monoxide	CO	10	0.05		
Nitrogen	N ₂	10	0.05	4	0.01
Ethane	C ₂ H ₆	1	0.005	2	0.01
Methanol	CH ₃ OH	0.1	0.005		
Carbon dioxide	CO ₂	1	0.005	1	0.005
Propane	C ₃ H ₈	1	0.005	0.1	0.005
i-Butane	i-C ₄ H ₁₀	1	0.005	0.1	0.005
n-Butane	n-C ₄ H ₁₀	1	0.005	0.1	0.005
i-Pentane	i-C ₅ H ₁₂	0.5	0.005	0.1	0.005
n-Pentane	n-C ₅ H ₁₂	0.5	0.005	0.1	0.005
n-Hexane	n-C ₆ H ₁₄			0.1	0.005
n-Heptane	n-C ₇ H ₁₆			0.1	0.005
HHV			0.3% relative		

Table 4 shows data from a Prima PRO factory test on a gravimetric cylinder containing 20 components, including inorganics and hydrocarbons from C₁ to C₆. The test was carried out over 18 hours and the measured standard deviations for all 20 components were lower than the specified standard deviations over 8 hours.

As well as the composition data, Prima PRO calculated the Compressibility, Ideal Lower Heating Value and Actual Lower Heating Value for the stream. Figure 6 shows these three parameters over the 18 hours of the test; as with the component concentrations, the measured LHV 18 hour standard deviations are well below the specified 8 hour standard deviations.

Table 4 Factory test on gravimetric cylinder containing 20 inorganic & hydrocarbon compounds, analyzed over 18 hours

	H ₂	CH ₄	C ₂ H ₂	N ₂	C ₂ H ₄	C ₂ H ₆	O ₂	H ₂ S (ppm)	Argon	C ₃ H ₆
Average	30.10	36.10	0.02	10.00	2.98	9.98	2.00	503.9	0.50	2.00
Cylinder Certificate	30	36.15	0.02	10	3	10	2	500	0.5	2
Actual Standard Deviation (18 hours)	0.023	0.015	0.0012	0.004	0.005	0.007	0.002	2.8	0.0005	0.002
Specified Standard Deviation (8 hours)	0.03	0.04	0.005	0.02	0.01	0.01	0.005	5	0.002	0.005
	CO ₂	C ₃ H ₈	1, 3 Butadiene	Butenes	iso Butane	n Butane	Pentenes	iso Pentane	n Pentane	n Hexane
Average	0.50	2.99	0.50	0.50	0.50	0.50	0.10	0.30	0.30	0.10
Cylinder Certificate	0.5	3	0.5	0.5	0.5	0.5	0.1	0.3	0.3	0.1
Actual Standard Deviation (18 hours)	0.003	0.002	0.0004	0.0008	0.0008	0.002	0.0004	0.001	0.002	0.001
Specified Standard Deviation (8 hours)	0.005	0.005	0.005	0.005	0.005	0.005	0.005	0.005	0.005	0.005

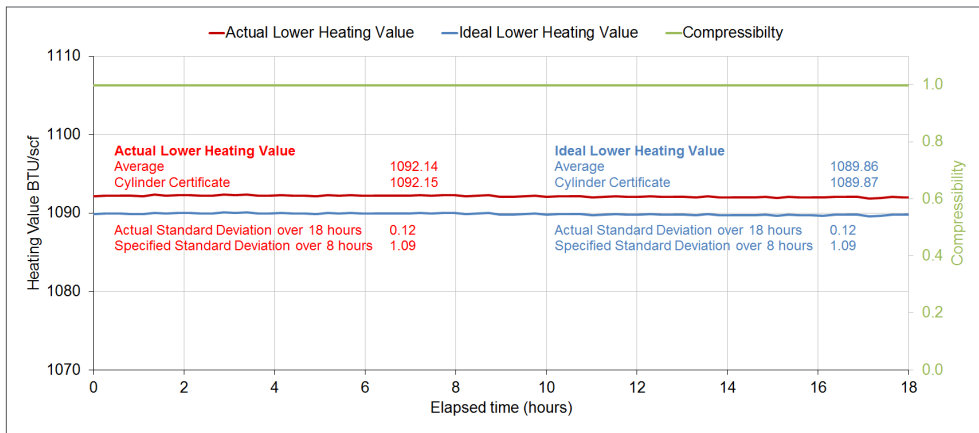


Figure 6 Heating values obtained during factory test on gravimetric cylinder containing 20 inorganic and hydrocarbon compounds, analyzed over 18 hours

Figure 7 shows 24 hours' process data from a flare gas stream. Prima PRO shows a spike in the ethylene concentration up to 5.8% after 3 hours, then a few small spikes of i-butane up to 5.5% after 12 hours, and finally a larger spike of i-butane up to 16.6% at around 23 hours. Identifying these individual components greatly assists the root cause analysis to locate the source of these emissions.

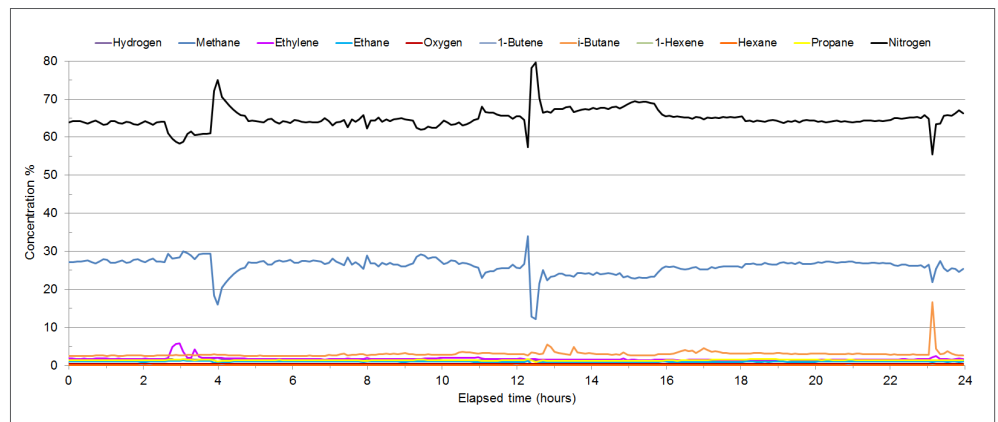


Figure 7 24-hour trend display of flare gas stream

Analysis of total sulfur

In general, regulators are interested in values for hydrogen sulfide (H_2S) and Total Reduced Sulfur (TRS). However there is some divergence on what constitutes TRS. In some cases it is defined as H_2S together with carbonyl sulfide (COS) and carbon disulfide (CS_2). It can also be defined as a mixture of compounds which contain a sulfur component in the reduced form, most commonly H_2S , methanethiol (methyl mercaptan, CH_3SH), dimethyl sulfide (DMS, $(CH_3)_2S$) and dimethyl disulfide (DMDS, $CH_3S_2CH_3$).

Mass spectrometers can measure a number of sulfur compounds down to ppm levels; typical performance figures are shown in Table 5. Analysis time is less than 30 seconds, including stream switching time and standard deviations are measured over 8 hours.

Table 5 Typical Prima PRO performance specification for sulfur compounds

Component	Typical Composition %mol	Precision of analysis by Prima PRO (single standard deviation) \leq
Hydrogen Sulfide	3 ppm	0.5 ppm
Methyl Mercaptan	10 ppm	0.5 ppm
Ethyl Mercaptan	10 ppm	0.5 ppm
n-Propyl Mercaptan	10 ppm	0.5 ppm
n-Butyl Mercaptan	10 ppm	0.5 ppm

Figure 8 shows 30 consecutive measurements on a test blend of three mercaptans at ppm levels. The standard deviations for the three compounds are shown in Table 6.

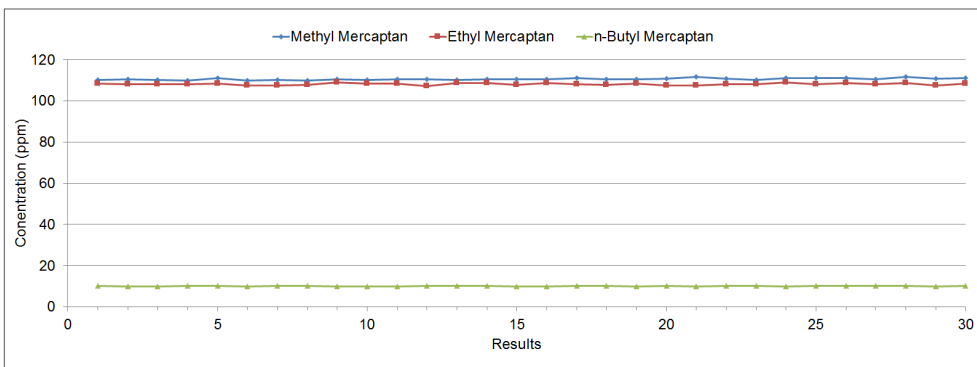


Figure 8 Test blend of three mercaptans

Table 6 Precision data for three typical sulfur compounds

	Methyl chloride	Vinyl chloride	Ethyl chloride	Ethylene dichloride
Mean	3.51	3.53	4.63	0.61
Standard deviation	0.04	0.06	0.16	0.05

Figure 9 demonstrates Prima PRO's excellent linearity for sulfur compounds. Eighteen hydrogen sulfide standards were generated by diluting a calibration cylinder using mass flow control; concentrations ranged from 4.85% down to 800 ppm. Part-per-million calibration standards can therefore be used with confidence to calibrate Prima PRO for sulfur compounds.

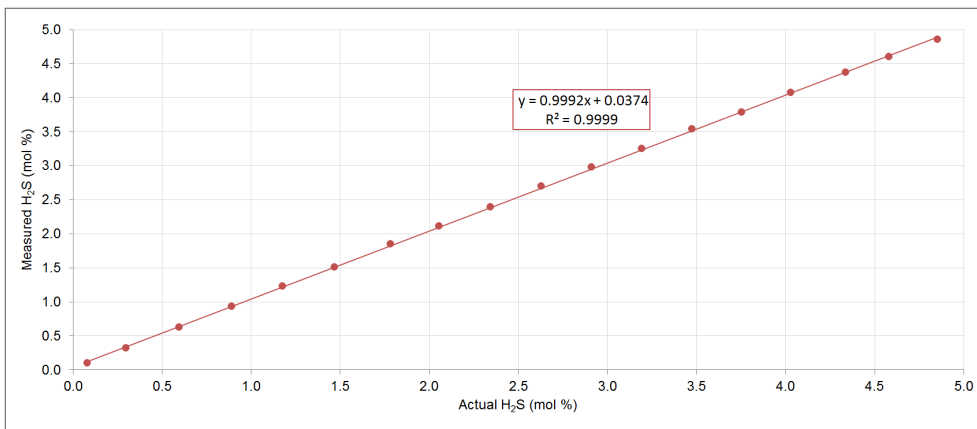


Figure 9 Linearity plot for H_2S from 800 ppm to 4.85%

GasWorks software can generate a figure for Total Sulfur using its Derived Value capability to sum the individual component concentrations. However it should be stressed that the total sulfur value only represents the sum of the sulfur compounds that the MS has analyzed—any unknown or unidentified compounds will not be reported.

For a true Total Sulfur reading we recommend the Thermo Scientific™ SOLA II™ Flare System which provides a solution for continuous and accurate determination of total sulfur in flare gas streams². SOLA II Flare uses PUVF (pulsed ultra-violet fluorescence) spectrometry to determine total sulfur. All organically bound sulfur is first converted to sulfur dioxide (SO₂) by sample combustion. Irradiation of SO₂ with ultraviolet light at a specific wavelength forms an excited form of SO₂. The excited SO₂ relaxes to its ground state by the emission of light or fluorescence. The intensity of the emitted light is directly proportional to the SO₂ concentration and thus the flare stack's total sulfur concentration.

Summary

Prima PRO Process MS provides fast on-line accurate analysis of flare gas composition. The inherent power of magnetic sector mass spectrometry, combined with the flexibility of GasWorks software, enables one Prima PRO to monitor not only flare gas streams, but also multiple process streams. Root cause fault analysis is therefore greatly facilitated, by comparing the detailed composition data from the flare gas stream with that of the various process streams.

As well as complete compositional analysis, Prima PRO provides accurate fuel gas properties, including LHV, HHV, density, specific gravity, Wobbe Index, stoichiometric air requirement and CARI. This ensures waste gases are burnt to complete combustion and unburned pollutants are not emitted from the flare.

Reference

1. EPA 40 CFR § 63.670—Requirements for Flare Control Devices, US Government Publishing Offices.
2. Continuous flare stack emission monitoring using the Thermo Scientific SOLA II Flare System Thermo Fisher Scientific Application Note, 2014.

USA

27 Forge Parkway
Franklin, MA 02038
Ph: (800) 437-7979
Fax: (713) 272-2273
orders.process.us@thermofisher.com

India

C/327, TTC Industrial Area
MIDC Pawane
New Mumbai 400 705, India
Ph: +91 22 4157 8800
india@thermofisher.com

China

+Units 702-715, 7th Floor
Tower West, Yonghe
Beijing, China 100007
Ph: +86 10 84193588
info.eid.china@thermofisher.com

Europe

Ion Path, Road Three,
Winsford, Cheshire CW73GA UK
Ph: +44 1606 548700
Fax: +44 1606 548711
sales.epm.uk@thermofisher.com

Find out more at thermofisher.com/flaregases

ThermoFisher
S C I E N T I F I C