

Ethylene Oxide and Propylene Oxide Analysis Using the Agilent 990 Micro GC

Authors

Fang Yue and Jiang Fei
Agilent Technologies
(Shanghai) Co. Ltd.

Abstract

This application note presents a method for the analysis of ethylene oxide (EO) and propylene oxide (PO) in their synthesis using an Agilent 990 Micro GC. Raw materials and products of interest (hydrogen [H₂], oxygen [O₂], nitrogen [N₂], carbon monoxide [CO], carbon dioxide [CO₂], ethylene, EO, propylene, and PO) before and after the reaction can all be separated and quantified on the Agilent J&W CP-Molsieve 5Å and Agilent J&W PoraPLOT U channels with excellent repeatability and a short analytical cycle time (less than 4 minutes). The lower detection limit (LDL) for EO and PO can be as low as parts per million (ppm).

Introduction

1,2-Epoxy compounds, abbreviated as epoxides, are a class of compounds with a ternary cyclic ether structure. The main epoxides in industry are EO and PO. Both are important chemical intermediates in the production of various solvents, surfactants, and polymer precursors, and they have a wide range of applications in the chemical, pharmaceutical, construction, pesticide, and other sectors.¹⁻³

The traditional synthesis of epoxides is carried out under high temperature and pressure, and the oxygen for epoxidation comes from the oxygen supply in the feed. In this situation, ethylene is prone to peroxidation, generating constant amounts of CO₂ and CO.¹⁻³ Thus, both raw materials and products are mostly constant components, and accurate analysis does not require high chromatographic performance.

However, in the electrocatalytic synthesis of epoxides, such as halogen-mediated indirect oxidation processes, the oxygen for epoxidation comes from water. The amount of free oxygen in the system, which is a byproduct of side reactions, is much less than that in traditional one. Furthermore, due to its mild reaction conditions, the peroxidation reaction of ethylene is significantly inhibited, and the yields of CO and CO₂ byproducts are much lower.^{1,3} Currently, this direction is still in the laboratory research stage. The overall reaction scale is small, with low current and yield, which puts high demands on the analytical equipment's detection limit and sample consumption.¹⁻³ At the same time, exploring reaction mechanisms efficiently and batch screening of catalysts requires rapid online analysis.

The Agilent 990 Micro GC can perform quick and accurate analysis of ppm-level EO, PO, and other related components from both traditional and/or electrocatalytic reaction systems. Its low consumption of power and carrier gas make it an ideal tool for online analysis, bringing value to both industrial production and laboratory research. Automatic sequential analysis of multiple samples and calibrations can also be achieved on one machine to accelerate batch screening by adding stream selection valves.

In this application note, two analytical channels of the 990 Micro GC were used, namely the 3+10 m Agilent J&W CP-Molsieve 5Å and the 1+10 m Agilent J&W PoraPLOT U. The 3+10 m Molsieve 5Å enabled the separation of permanent gases, while the 1+10 m PoraPLOT U enabled the separation of CO₂, ethylene, and EO, or CO₂, propylene, and PO under corresponding conditions.⁴ Each run took less than 4 minutes to complete, with excellent repeatability for multiple analyses.

Experimental

The 990 Micro GC was equipped with a 3+10 m Molsieve 5Å backflush channel (retention time stability configuration), and a 1+10 m PPU backflush channel. Table 1 shows the experimental conditions for the analysis. Standard gases were purchased from Air Liquide, Inc. (Shanghai, China). Table 2 shows the composition of the standard gases.

Table 1. Experimental conditions for EO and PO synthesis analysis using the 990 Micro GC.

| Application Scope | EO and PO Synthesis | EO Synthesis | PO Synthesis |
|----------------------|-------------------------------|------------------------------|------------------------------|
| Channel Type | 3+10 m Molsieve 5Å, backflush | 1+10 m PoraPLOT U, backflush | 1+10 m PoraPLOT U, backflush |
| Carrier Gas | Helium | Helium | Helium |
| Column Pressure | 150 kPa | 150 kPa | 125 kPa |
| Injector Temperature | 65 °C | 70 °C | 70 °C |
| Column Temperature | 65 °C | 70 °C | 110 °C |
| Injection Time | 40 ms | 40 ms | 40 ms |
| Backflush Time | 15.8 s | 14.0 s | 17.8 s |
| Sample Time | 60 s | 60 s | 60 s |

Table 2. Composition of the standard gases.

| Standard Gas 1 | | Standard Gas 2 | |
|-----------------|---------------|-----------------|---------------|
| Component | Concentration | Component | Concentration |
| N ₂ | Balance | N ₂ | Balance |
| Ethylene | 75.0% | Propylene | 74.7% |
| H ₂ | 390 ppm | H ₂ | 410 ppm |
| O ₂ | 194 ppm | O ₂ | 200 ppm |
| EO | 90.0 ppm | PO | 114 ppm |
| CO | 50.9 ppm | CO | 52.0 ppm |
| CO ₂ | 49.8 ppm | CO ₂ | 47.5 ppm |

Results and discussion

The Molsieve 5Å is used for the separation of permanent gases. To adapt to different sample compositions, especially for high-concentration CO₂ analysis in traditional synthesis processes, a 3-meter-long precolumn is preferred for the Molsieve 5Å channel to ensure that CO₂ will not enter the analysis column. Based on the actual sample system, other Molsieve 5Å configurations may also be selected. Helium is applied as the carrier gas to elicit a higher response in components other than H₂, and the actual content of H₂ in the sample is unlikely to exceed 10%, so there is no need to worry about peak inversion brought on by the abnormal thermal conductivity of the H₂/He mixture.⁵ However, due to the close thermal conductivity between He and H₂, analysis of H₂ component with a concentration lower than 50 ppm is unlikely.

The PoraPLOT U channel is used for the separation of permanent gas, CO₂, olefin raw materials, and the target product (EO or PO). A straight configuration can also be chosen. The use of backflush configuration is recommended here to allow any possible higher-boiling components in these complex practical systems to be vented out from the column, keeping analysis time short. Since the synthesis of EO and PO is generally independent of each other, and the retention time (RT) of EO and PO on the PoraPLOT U under the same

conditions differs greatly, different analysis conditions were applied for EO and PO to shorten analysis time, enhance sensitivity, and improve separation. Due to the significantly stronger retention of epoxides on the PoraPLOT U, the overall principle for adjusting analysis conditions is to minimize the RT of epoxides while ensuring the effective separation of other components. The separation of CO₂ and ethylene in the EO system and the separation of permanent gas and CO₂ in the PO system are critical for the respective analyses.

The analysis results of the EO synthesis were as follows. Figure 1A shows the chromatogram of Standard Gas 1 on the 3+10 m Molsieve 5Å backflush channel. H₂, O₂, N₂, and CO were well resolved. On the 1+10 m PoraPLOT U backflush channel, permanent gas composite, CO₂, ethylene, and EO were separated as shown in Figure 1B. EO had a good peak shape.

The analysis results of the PO synthesis were as follows. Figure 2A shows the chromatogram of Standard Gas 2 on the 3+10 m Molsieve 5Å backflush channel. H₂, O₂, N₂, and CO were well resolved, as expected. On the 1+10 m PoraPLOT U backflush channel, permanent gas, CO₂, ethylene, and PO were separated as shown in Figure 2B. PO also had a good peak shape.

The maximum analytical cycle time was less than 4 minutes for both the EO and PO methods.

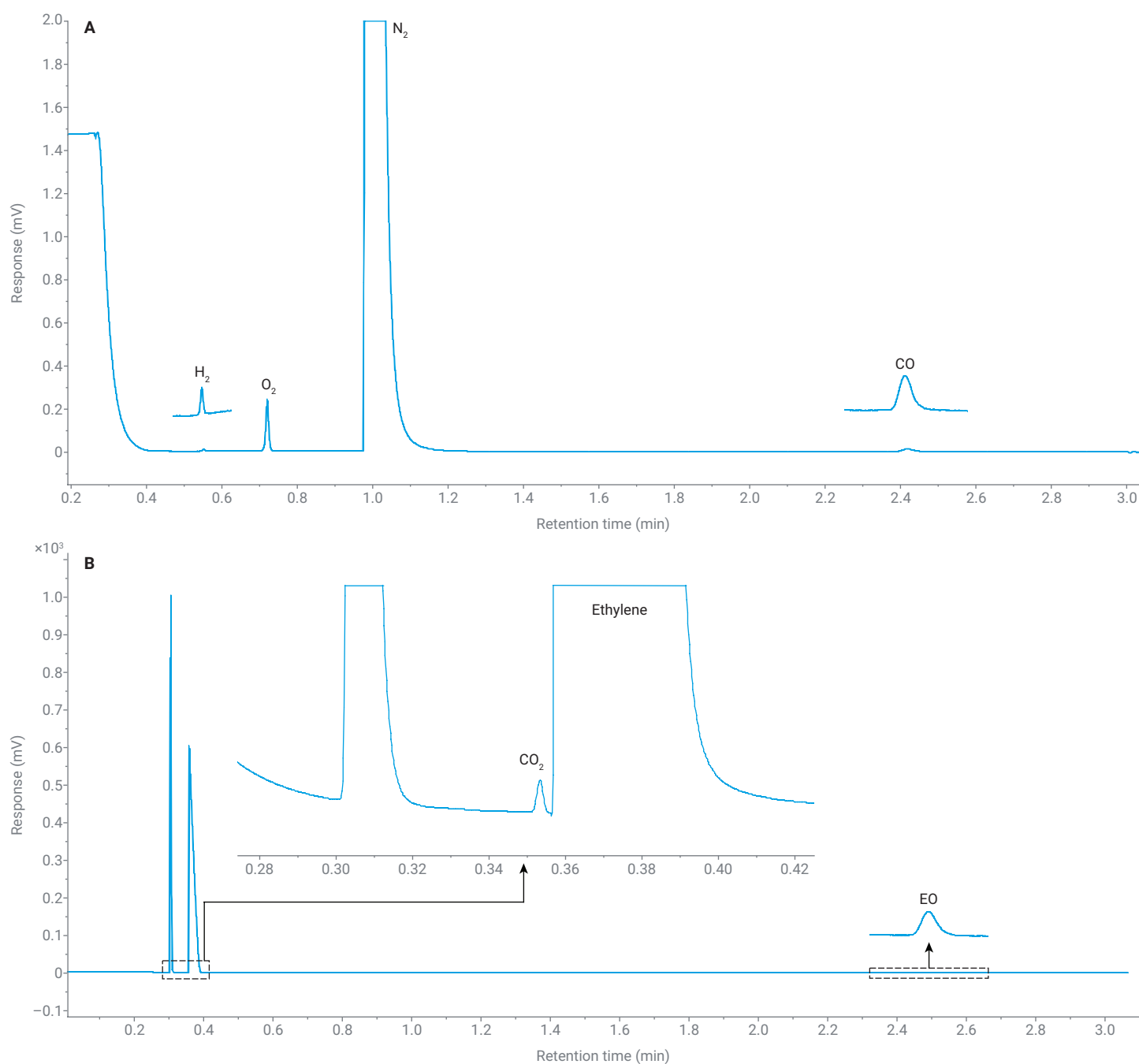


Figure 1. (A) Chromatogram of the Standard Gas 1 (EO synthesis) on the 3+10 m Agilent J&W CP-Molsieve 5Å backflush channel. (B) Chromatogram of the Standard Gas 1 (EO synthesis) on the 1+10 m Agilent J&W PoraPLOT U backflush channel.

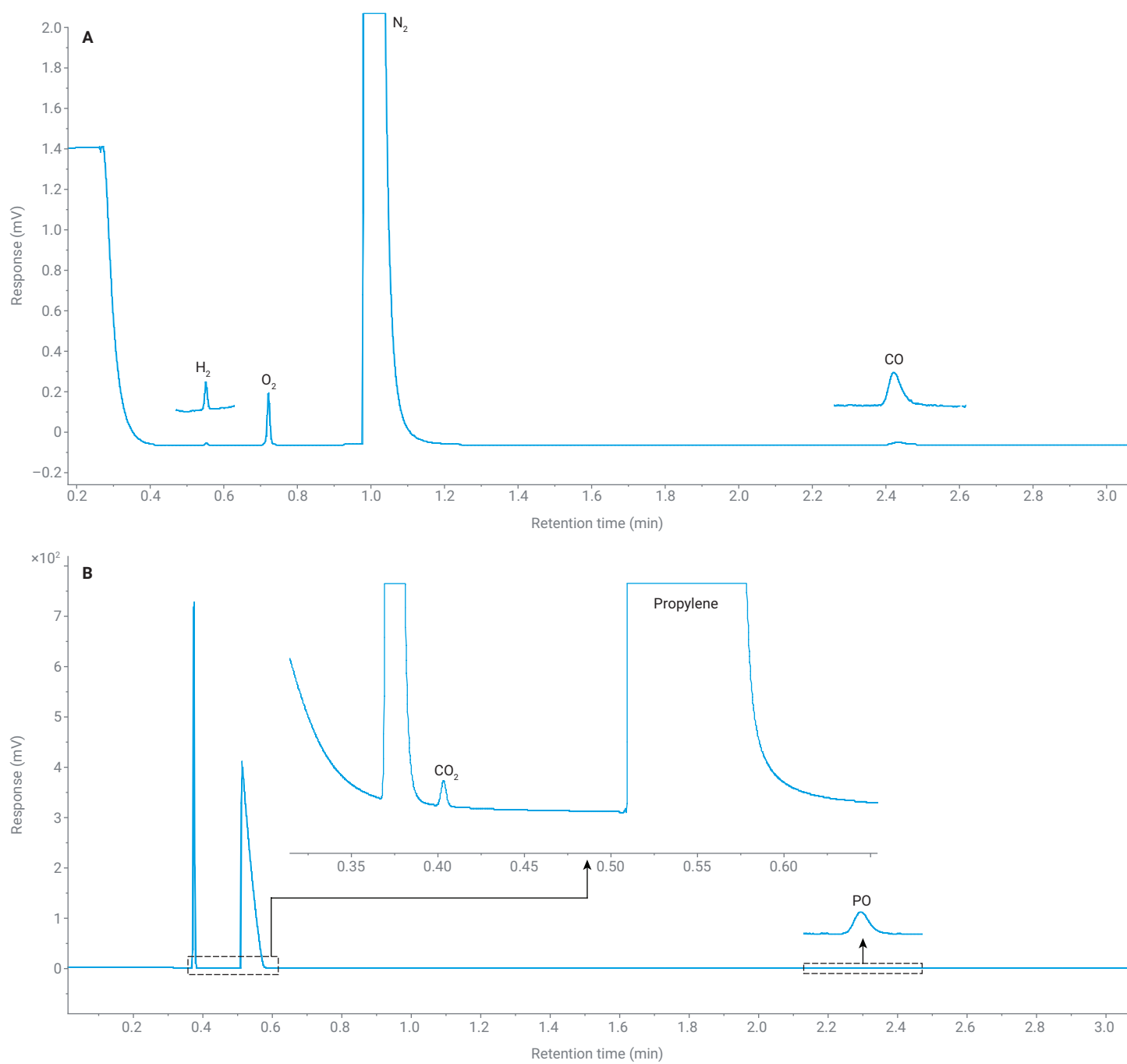


Figure 2. (A) Chromatogram of the Standard Gas 2 (PO synthesis) on the 3+10 m Agilent J&W CP-Molsieve 5Å backflush channel. (B) Chromatogram of the Standard Gas 2 (PO synthesis) on the 1+10 m Agilent J&W PoraPLOT U backflush channel.

Table 3 and Table 4 show the performance for a sequence of 10 runs of these two standard gases. The RT and area relative standard deviations (%RSDs) for these components are less than 0.2% and 2.5%, respectively. These low %RSD values demonstrate the excellent performance of the 990 Micro GC and its ability to provide quantitative results with a high level of confidence, repeatability, and quality.

Table 5 shows the estimated LDL of related components on the 3+10 m Molsieve 5Å backflush and 1+10 m PoraPLOT U backflush based on the calculation of current results or historical data that was obtained under similar experimental conditions described in Table 1.

Table 3. Retention time, area, and repeatability (%RSD) of these measures for 10 runs of the Standard Gas 1 (EO synthesis).

| Component | Concentration | RT (min) | RT RSD (%) | Area (mV × s) | Area RSD (%) | Analysis Channel |
|-----------------|----------------|----------|------------|---------------|--------------|----------------------------------------------|
| H ₂ | 390 ppm | 0.552 | 0.011 | 0.005 | 2.428 | 3+10 m Agilent J&W CP-Molsieve 5Å, backflush |
| O ₂ | 194 ppm | 0.720 | 0.005 | 0.125 | 0.526 | 3+10 m Agilent J&W CP-Molsieve 5Å, backflush |
| N ₂ | Balance (~25%) | 0.985 | 0.012 | 156.582 | 0.017 | 3+10 m Agilent J&W CP-Molsieve 5Å, backflush |
| CO | 50.9 ppm | 2.419 | 0.040 | 0.032 | 2.137 | 3+10 m Agilent J&W CP-Molsieve 5Å, backflush |
| CO ₂ | 49.8 ppm | 0.354 | 0.040 | 0.034 | 0.141 | 1+10 m Agilent J&W PoraPLOT U, backflush |
| Ethylene | 75.0% | 0.359 | 0.000 | 516.881 | 0.093 | 1+10 m Agilent J&W PoraPLOT U, backflush |
| EO | 90.0 ppm | 2.493 | 0.037 | 0.067 | 1.691 | 1+10 m Agilent J&W PoraPLOT U, backflush |

Table 4. Retention time, area, and repeatability (%RSD) of these measures for 10 runs of the Standard Gas 2 (PO synthesis).

| Component | Concentration | RT (min) | RT RSD (%) | Area (mV × s) | Area RSD (%) | Analysis Channel |
|-----------------|----------------|----------|------------|---------------|--------------|----------------------------------------------|
| H ₂ | 410 ppm | 0.553 | 0.005 | 0.005 | 2.540 | 3+10 m Agilent J&W CP-Molsieve 5Å, backflush |
| O ₂ | 200 ppm | 0.722 | 0.003 | 0.132 | 0.174 | 3+10 m Agilent J&W CP-Molsieve 5Å, backflush |
| N ₂ | Balance (~25%) | 0.986 | 0.006 | 163.553 | 0.018 | 3+10 m Agilent J&W CP-Molsieve 5Å, backflush |
| CO | 52.0 ppm | 2.433 | 0.047 | 0.032 | 2.556 | 3+10 m Agilent J&W CP-Molsieve 5Å, backflush |
| CO ₂ | 47.5 ppm | 0.404 | 0.004 | 0.037 | 0.102 | 1+10 m Agilent J&W PoraPLOT U, backflush |
| Propylene | 74.7% | 0.513 | 0.000 | 717.113 | 0.041 | 1+10 m Agilent J&W PoraPLOT U, backflush |
| PO | 114 ppm | 2.309 | 0.041 | 0.078 | 0.940 | 1+10 m Agilent J&W PoraPLOT U, backflush |

Table 5. Lower detection limit (LDL) of this configuration.

| Component | LDL (ppm) | Analysis Channel |
|-----------------|-----------|----------------------------------------------|
| H ₂ | 50 | 3+10 m Agilent J&W CP-Molsieve 5Å, backflush |
| O ₂ | 10 | 3+10 m Agilent J&W CP-Molsieve 5Å, backflush |
| N ₂ | 10 | 3+10 m Agilent J&W CP-Molsieve 5Å, backflush |
| CO | 5 | 3+10 m Agilent J&W CP-Molsieve 5Å, backflush |
| CO ₂ | 1 | 1+10 m Agilent J&W PoraPLOT U, backflush |
| Ethylene | 2 | 1+10 m Agilent J&W PoraPLOT U, backflush |
| EO | 5 | 1+10 m Agilent J&W PoraPLOT U, backflush |
| Propylene | 2 | 1+10 m Agilent J&W PoraPLOT U, backflush |
| PO | 5 | 1+10 m Agilent J&W PoraPLOT U, backflush |

Conclusion

This application note demonstrates the chemical performance of Agilent 990 Micro GC for the analysis of ethylene oxide (EO) and propylene oxide (PO) synthesis. The 3+10 m Agilent J&W CP-Molsieve 5Å and the 1+10 m Agilent J&W PoraPLOT U channel on the 990 Micro GC can analyze H₂, O₂, N₂, CO, CO₂, ethylene, EO, propylene, and PO with excellent repeatability, a short analytical cycle time, and sufficient LDL at ppm levels. The 990 Micro GC is well suited for the fast and precise analysis of EO, PO, and their synthesis-related components.

References

1. Wan Ru, L. Chloride-Mediated Selective Electrosynthesis of Ethylene and Propylene Oxides at High Current Density. *Science* **2020**, 368, 1228–1233.
2. Yixuan, S. Research Progress on Electrochemical Synthesis of Ethylene Oxide and Propylene Oxide. *Petrochemical Technology* **2023**, 52(11), 1596–1601.
3. Hanyu, W.; *et al.* Boosting Electrocatalytic Ethylene Epoxidation by Single Atom Modulation. *Angew. Chem. Int. Ed.* **2024**, 63, e202402950.
4. Jie, Z. Fast Analysis of Natural Gas using the Agilent 990 Micro GC Natural Gas Analyzer, *Agilent Technologies application note*, publication number 5994-1040EN, **2019**.
5. Toonen, A.; van Loon, R. Hydrogen Detection with a TCD using Mixed Carrier Gas on the Agilent Micro GC, *Agilent Technologies application note*, publication number 5991-3199EN, **2013**.