

Dynamic gas-solid surface interactions for dry reforming of methane via in-situ DRIFTS analysis

Authors

Md Monir Hossain, Ruigang Wang,
Michigan State University

In this note: Monitoring surface species on a nanorod catalyst for dry reforming of methane

Dry reforming of methane (DRM) is a key process for converting CH_4 and CO_2 into valuable syngas (CO and H_2).¹⁻³ Understanding gas-solid surface interactions during DRM is essential for the rational design of catalysts.

Here, in situ Diffuse Reflectance Infrared Fourier Transform Spectroscopy (DRIFTS) using the Harrick Praying Mantis™ accessory and a high-temperature reaction chamber (HVC-DRM-5) are employed to track surface species on a 1.5 wt% Ni–0.5 wt% Ru/ CeO_2 nanorod (NR) catalyst over a temperature range of 50°C to 450°C.

Experimental

Catalyst preparation

The 1.5 wt% Ni–0.5 wt% Ru/ CeO_2 NR catalyst was synthesized via a one-pot hydrothermal method, ensuring uniform dispersion of Ni–Ru on the CeO_2 NR supports.

Instrumentation

In situ DRIFTS was performed with the Harrick Praying Mantis diffuse reflection accessory along with a high-temperature reaction chamber, integrated into a Thermo Scientific™ Nicolet™ iS50 FTIR Spectrometer system (Fig. 1), providing precise temperature and gas flow control for real-time monitoring of surface gas-solid reactions.

Reaction conditions

A CH_4/CO_2 gas mixture (29:71 $\approx$ 0.4 ratio) was fed at a total flow rate of 100 sccm. The temperature was ramped from 50°C to 450°C at 10°C/min, with spectra collected every 50°C after stabilizing the gas flow and temperature for 15 min.

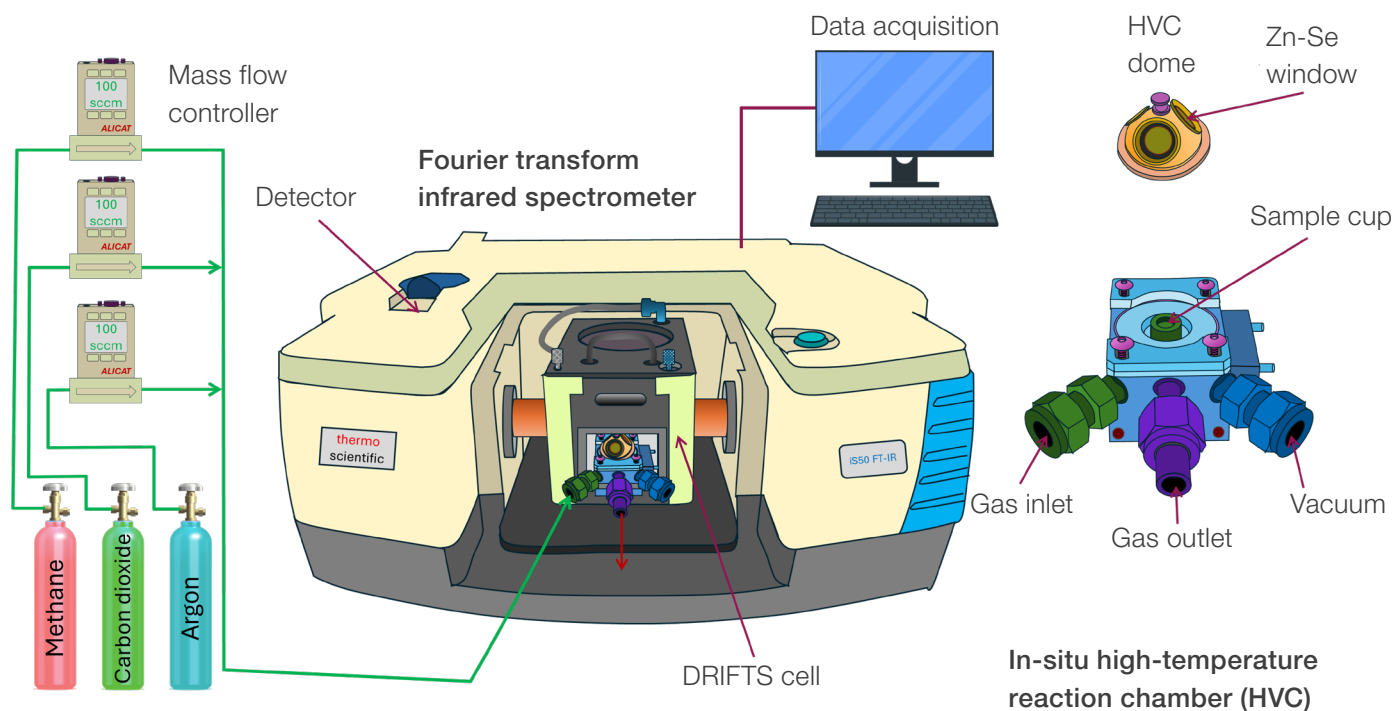


Figure 1. Experimental setup for in-situ DRIFTS using a Harrick Praying Mantis Cell and a high-temperature reaction chamber.

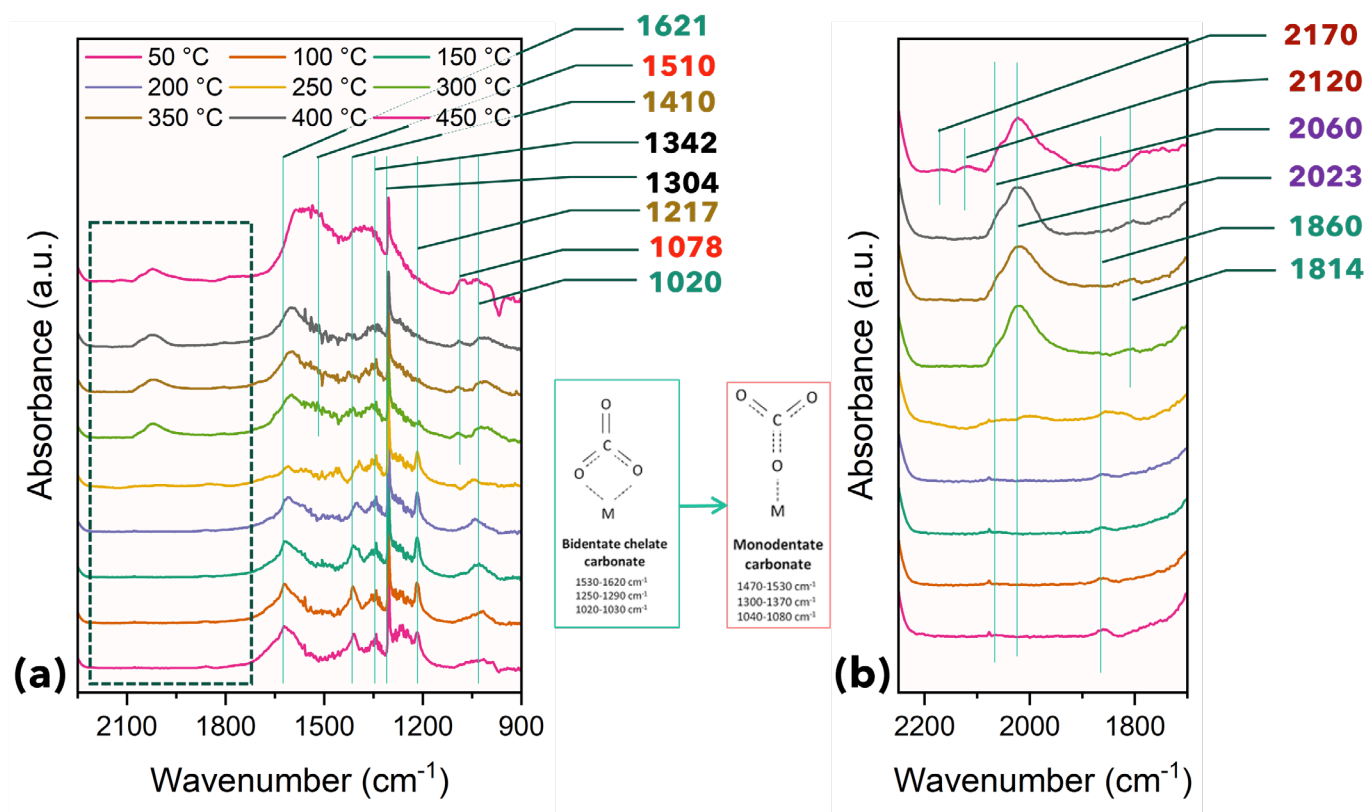


Figure 2. In situ DRIFTS spectra for DRM at varying temperatures ($\text{CH}_4/\text{CO}_2 = 29/71$ sccm ≈ 0.4 ratio; total flow rate: 100 sccm).

Conclusion

The in situ DRIFTS investigation with the Harrick Praying Mantis setup revealed the evolution and transformation of surface species during DRM on Ni–Ru/CeO₂ NRs. The findings demonstrate the power of in situ DRIFTS's ability to elucidate catalytic mechanisms and highlight the promising performance of this catalyst for DRM.

Acknowledgement

This project is supported by the U.S. National Science Foundation (2427238 and 2427213).

References

1. Ahasan, M. R.; Hossain, M. M.; Wang, R. Dielectric Barrier Discharge Reactors for Plasma-Assisted CO₂ and CH₄ Conversion: A Comprehensive Review of Reactor Design, Performance, and Future Prospects. *Energy Technology* 2024, 13 (4), 2401177.
2. Monir Hossain, M.; Robayet Ahasan, M.; Wang, R. Influence of Catalyst Shape on Plasma-Assisted Dry Reforming of Methane: A Comparative Study of Ni–CeO₂ Nano-Cubes and Nano-Octahedra. *Chemical Engineering Journal* 2024, 496, 154193.
3. Hossain, M. M.; Ahasan, M. R.; Wang, R. Exploring the Versatile Roles of Microwave Energy in Catalyst Synthesis, Treatment, and Catalytic Performance. *Fuel* 2025, 400, 135684.
4. Vakili, R.; Gholami, R.; Stere, C. E.; Chansai, S.; Chen, H.; Holmes, S. M.; Jiao, Y.; Hardacre, C.; Fan, X. Plasma-Assisted Catalytic Dry Reforming of Methane (DRM) over Metal–Organic Frameworks (MOFs)-Based Catalysts. *Appl Catal B* 2020, 260, 118195.
5. Cárdenas-Arenas, A.; Quindimil, A.; Davó-Quinonero, A.; Bailón-García, E.; Lozano-Castelló, D.; De-La-Torre, U.; Pereda-Ayo, B.; González-Marcos, J. A.; González-Velasco, J. R.; Bueno-López, A. Isotopic and in Situ DRIFTS Study of the CO₂ Methanation Mechanism Using Ni/CeO₂ and Ni/Al₂O₃ Catalysts. *Appl Catal B* 2020, 265, 118538.
6. Li, Y.; Qin, T.; Wei, Y.; Xiong, J.; Zhang, P.; Lai, K.; Chi, H.; Liu, X.; Chen, L.; Yu, X.; Zhao, Z.; Li, L.; Liu, J. A Single Site Ruthenium Catalyst for Robust Soot Oxidation without Platinum or Palladium. *Nature Communications* 2023 14:1 2023, 14 (1), 1–11.

 Learn more at thermofisher.com/is50

thermo scientific

For research use only. Not for use in diagnostic procedures. For current certifications, visit thermofisher.com/certifications

© 2026 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. Harrick Praying Mantis is a registered trademark of Specac Ltd.

MCS-AN1911-EN 0826