ChemiSorb HTP

High Thoughput Chemisorption Analyzer

- High throughput with six analysis stations
- Up to six chemisorption analyses can begin simultaneously
- Independent furnaces for each analysis station features a user defined temperature range from 10 °C above ambient to 700 °C.
- Quartz sample reactor with flow-through design, available for various size pellets, core and powders
- Fully automated analysis
- Long unattended analysis times for high resolution adsorption isotherms
- Analysis ports operate concurrently or in parallel
- Up to 12 different gases can be attached simultaneously
- Windows® driven software

Optimum design and efficient utilization of catalysts require a thorough understanding of the surface structure and surface chemistry of the catalytic material. Chemical adsorption analyses can provide much of the information needed to evaluate catalyst materials in the design and production phases, as well as after a period of use.

The ChemiSorb HTP is a fully-automated high-throughput chemisorption analyzer that determines the percent metal dispersion, active metal surface area, size of active particles, and surface acidity of catalyst materials. A major advantage of this instrument is its six analysis stations. Multiple analyses can be run simultaneously or in parallel on one instrument not only saving time for busy catalyst operations, but providing economy of lab space. The ChemiSorb HTP features a vacuum system and manifold with constantly monitored pressure transducers on each of its six analysis stations. Each port is also equipped with a furnace enabling independent control of sample temperature and ramping from 10 °C above ambient to 700 °C. The temperature is recorded with each equilibrium pressure to provide the highest quality isotherms possible. An equilibration option allows the user to specify different equilibration times. A mass flow controller is installed in each port to ensure accurate and reproducible flow through the sample. Up to twelve different gases can be attached to the ChemiSorb HTP simultaneously.

Degas/sample preparation is done in-situ prior to analysis. Samples may be added or removed from each station without



disturbing the treatment of other samples undergoing preparation or analysis. The instrument features long unattended analysis times and a high-throughput mode that allows the user to start multiple parallel analyses.

The ChemiSorb HTP analysis program operates in a Windows[®] environment. This makes operation of the analyzer easier and allows the user to run other applications while an automatic operation is in progress. The report system provided in the analysis program allows the user to manipulate and customize reports in a variety of ways. Zoom in on portions of the graphs or shift the axes to examine fine details. Graphs and data can be copied to the clipboard and pasted into other applications. Reports can be customized with a choice of fonts and a company logo added to the report header.



www.micromeritics.com

Micromeritics Instrument Corporation 4356 Communications Dr., Norcross, GA 30093, USA US Sales: 770.662.3633, International Sales: 770.662.3660 Customer Orders: 770.662.3636, FAX: 770.662.3696

Specifications

Analysis **Furnace System** Capacity: 6 sample stations The ChemiSorb HTP has six independent furnaces, one for each sample port. **Manifold Temperature** Seven platinum resistance detector (RTD) Type: Temperature Range: 10 °C above ambient to 700 °C Accuracy: ± 0.02 °C by keyboard entry ± 1.0 °C Accuracy: ± 0.1 °C per month Stability: Stability: ± 0.5 °C up to 50 °C; **Gas Flow Control** \pm 1.0 °C from 50 to 700 °C; Type: Six Mass Flow Controller, One per port Ramp Range: 1 to 20 °C/min up to 500 °C 0 to 200 sccm (nitrogen equivalent flow rate) Range: 1 to 10 °C from 500 to 700 °C Accuracy: \geq 5% full scale Cool down: From 700 to 10 °C above **Gas Dosing** ambient in less than 60 minutes 0.05 cm3 STP, normal doses Accuracy Temperature increment: 1 °C 0.01 cm³ STP, low-pressure doses **Electrical Pressure Measurement** 100/115/230 VAC Voltage: Analysis Manifold Transducers: Frequency: 50 or 60 Hz Range: 0 to 950 mmHg Power: 1500 VA. maximum Resolution: 0.001 mmHg, 1000-mmHg transducer (exclusive of vacuum fore pump which is 0.00001 mmHg, 10-mmHg transducer powered separately) Accuracy: Within 0.15% of reading, 1000- and Environment 10-mmHg transducers (Includes Temperature: 10 to 35 °C (50 to 95 °F), operating; nonlinearity, hysteresis and non--10 to 55 °C (14 to 131 °F), storing or repeatability; transducer manufacturer's shipping specifications.) Humidity: Up to 90% non-condensing Sample Port Transducers: (for instrument) Range: 0 to 1000 mmHg 20 to 80% (for computer system) Resolution: 0.001 mmHg, 1000-mmHg transducer **Physical** 0.00001 mmHg, 10-mmHg transducer Width: 159 cm (63 in.) Within 0.5% of reading, 1000- and Accuracy: (20.2 in.) Depth: 51 cm 10-mmHg transducers (Includes non Height: 159 cm (62.5 in.) linearity, hysteresis and non repeatability; Weight: 215 kg (475 lb) transducer manufacturer's specifica-Computer tions.) Minimum requirements: Computer capable of running Windows® Vacuum Gauge: 7, Windows[®] XP Professional or Type: Pirani thermal vacuum gauge Windows Vista® Business or Ultimate 0.000375 mmHg to \geq 1 mmHg Range: operating system Resolution on screen: CD-ROM drive 0.0001 mmHg (0.1 microns) 512 megabytes of main memory 20-gigabyte hard drive Vacuum System SVGA monitor (1024 x 768 minimum Diaphragm pump: Two-stage, ultimate vacuum less than 3 mmHg resolution) High-vacuum pump: Less than 3.8×10^{-9} mmHg, measured by pump Ethernet port, capable of communicating manufacturer according to Pneurop Standard 5608 with a 10 base-T ethernet card

In keeping with a policy of ongoing product improvement, specifications are subject to change without notice



QUALITY MANAGEMENT SYSTEM CERTIFIED BY DNV Micromeritics Instrument Corporation 4356 Communications Dr., Norcross, GA 30093, USA US Sales: 770.662.3633, International Sales: 770.662.3660 Customer Orders: 770.662.3636, FAX: 770.662.3696

www.micromeritics.com