



ANATEL PAT700

TOTAL ORGANIC CARBON AND CONDUCTIVITY ANALYZER

Fully Compliant with Global Pharmacopeia Requirements : **USP, EP, JP**



CHARACTERIZED
by ingenuity.



PAT700 TOTAL ORGANIC CARBON & CONDUCTIVITY ANALYZER

LOW COST OF OWNERSHIP

TOC, conductivity and water temperature from just one analyzer

- Can be fully validated for TOC, Conductivity and Temperature to USP, EP and JP requirements

12 month service interval

- Auto-switching main and standby UV lamps
- UV Detect to ensure UV lamp is working correctly

No chemicals

- Designed specifically for pharmaceutical PW and WFI measurement, this analyzer uses just powerful UV light to oxidize the organics

No peristaltic pump

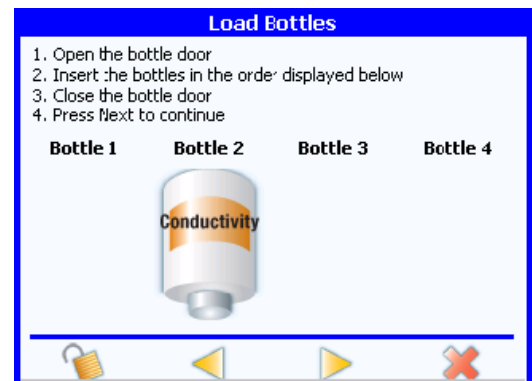
- No tubing and pump heads to replace
- PAT700 traps an aliquot for each analysis, so the measurement is stable and not affected by changes in sample pressure or flow rate

Single sensor to measure TIC and TC

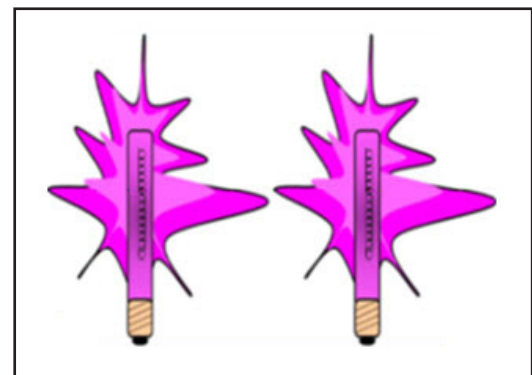
- $TC - TIC = TOC$
- Stable measurements for >12months

On-line and grab-sample analysis in one analyzer

- Built-in, 4-bottle grab-sample analyzer



Fully compliant conductivity meter



UV Monitoring with auto-switching
Main & Standby UV



IMPROVED COMPLIANCE

Fully ICH Q2 compliant

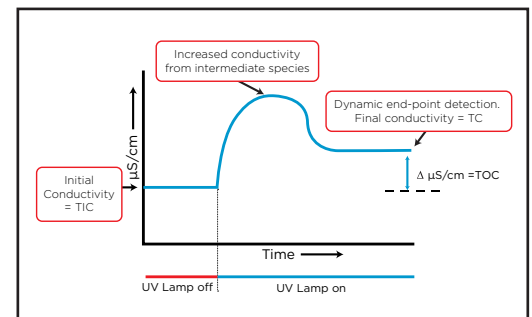
- Not affected by interference from ppm levels of TIC

Complete oxidation

- Fully compliant with EP 2.2.44 requirement for complete TOC oxidation through dynamic end-point detection technology

Root cause analysis support

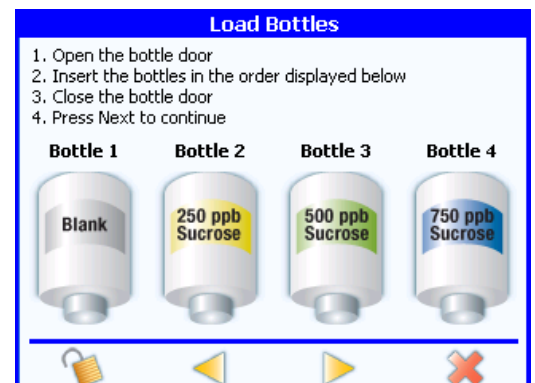
- Excursion capture feature allows a water sample to be captured to assist root cause analysis should a TOC excursion be detected
- Built-in grab-sample analyzer for analyzing samples from other points in the water loop



Complete TOC oxidation to EP2.2.44

21 CFR Part 11 Compliant

- Multi-level user access
- Windows credentials (Microsoft Active Directory)
- Secure .pdf file export via FTP over Ethernet
- No manual data entry – calibration standards use RFID tags to transfer lot number, expiry date and certified value directly to the PAT700
- Built-in electronic automated Calibration and System Suitability SOPs
- No manual calculations – automated Calibration and System Suitability Pass/Fail calculations



Built-in, automated SOPs

Specifications

TOC	Operating Range	0.5 to 2,000 ppb as Carbon
	Display Resolution	0.1 ppb
	Accuracy	±3 ppb or ±5%, whichever is greater
	Repeatability	±0.3 ppb or ±1%, whichever is greater
	Limit of Detection	0.5 ppb
	Maximum Input Conductivity	0.2 µS/cm for all waters, 1.0 µS/cm for all neutral waters, 5.0 µS/cm for water with CO ₂ as the sole conductive species
Conductivity	Conductivity	Range 0.05 to 150 µS/cm (@ 25°C)
	Display Resolution	0.01 µS/cm
	Conductivity Accuracy	±2% over full range (uncompensated)
	Available Conductivity Reporting Modes	Temperature compensated to 25°C, or uncompensated
	Available Resistivity Reporting Mode	Temperature compensated to 25°C only
	Resistivity	Resistivity Range 0.2 to 18 MΩ-cm (@ 25°C)
	Display Resolution	0.01 over full range
Temperature	Ambient Operating Range	10 to 40°C (50 to 104°F)
	Measurement Accuracy	±0.5°C
	Sample Water Range	1 to 95°C (34 to 203°F)
	Display Resolution	0.1 over full range
Physical Specs	UV Lamps	2, with UV Detect technology
	Interface/Display	Color touchscreen
	Maximum Altitude	4,000 m (13,125 ft)
	User I/O Wiring	Three, ¾-inch conduit openings or quick disconnect fittings
	Standards System	Onboard, Automated Standards Introduction System (OASIS)
	Dimensions	59.7 w X 22.9 d X 25.4 h cm (23.5 X 9 X 10 inches)
	Weight	13.6 kg (30 lbs)
	Sample Inlet Flow Rate Range	60 mL/min to 300 mL/min
	Sample Inlet Pressure Range	10 to 100 psi (69 to 690 kPa)
Communications	Analog output	3 x 4-20mA outputs, user configurable TOC, Conductivity (uncompensated) and Sample Temperature
	Digital output	4 x digital outputs, user configurable (for alarms, etc.)
	Digital input	2 x digital inputs (for remote control)
Compliance	Installation Category	II
	Pollution Degree	2, IEC 61010-1
	CE Compliance	EN 61010-1 and EN 61326
	Safety Rating	ETL, conforming to UL 61010-1 and CSA 22.2 No. 61010-1
	Enclosure Rating	Conduit version: IP56 Quick connect version: IP46
	Release tests,	USP <643>, USP <645>, JP 2.59, EP 2.2.44
New Features	CIP	Selectable mode for Clean-In-Place analysis
	Dual Stream option	Toggle or programmable stream switching
	Excursion sampling	Minimum flow rate to fill excursion bottle = 160 mL/min
	Rouge detection	Identifies oxidation cell contamination from rouging



© 2022 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo, and the Beckman Coulter product and service marks mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries. All other trademarks are property of their respective owners.

For Beckman Coulter's worldwide office locations and phone numbers, please visit Contact Us at beckman.com

22.06.5119.PCC