

# ANATEL PAT700 TOC ANALYZER PLATINUM AND GOLD



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



Facilities have different needs, refer to this comparison table to see which configuration best meets those needs.

The PAT700 is specifically designed to help demonstrate compliance to the pharmacopeial requirements for TOC and conductivity for Purified Water and Water For Injection.

ITEM NO.	DESCRIPTION
PAT700210	<b>PAT700 Platinum</b> PAT700200 - MULTIBOTTLE ANALYSIS TOC, US/JAPAN, QC (RFID)
PAT700215	<b>PAT700 Gold</b> PAT700200 - TOC, US/JAPAN, QC (NO RFID)

FEATURES	PAT700 PLATINUM	PAT700 GOLD
User Name and Password Controlled by Microsoft Active Directory	■	
Water sample capture in case of TOC Excursion	■	
Automatic RFID Recognition of Lot ID, calibration standards and System Suitability	■	
Manual Entry of Standards Information		■
21 CFR Part 11 Compliance Tools	■	■
Number of Users = 100	■	■
USP <643> Complete TOC Oxidation USP <645> Conductivity	■	■
12 month service interval: <ul style="list-style-type: none"> <li>Auto-switching main standby UV lamps</li> <li>UV Detect to ensure UV lamp is working correctly</li> </ul>	■	■
Built-in Pump and Heat Exchanger	■	■
Conduit Connections available	■	

# SPECIFICATIONS

---

<b>TOC</b>	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 <sub>2</sub> as the sole conductive species

---

<b>Conductivity</b>	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	

---

<b>Temperature</b>	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

---

<b>Physical Specs</b>	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)	

---

<b>Communications</b>	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)

---

<b>Compliance</b>	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

---

<b>New Features</b>	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

\* Not available on PAT700 Gold, non-RFID model



© 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](http://beckman.com).  
22.06.5118.PCC