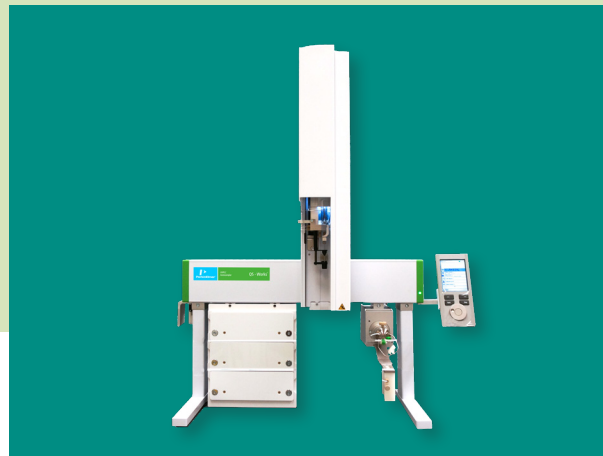


Liquid Chromatography/  
Mass Spectrometry

## QS-Works Autosampler



The QS-Works™ UHPLC Autosampler is part of PerkinElmer's QSight LC-MS/MS solution. This Autosampler provides enhanced performance, large sample capacity, high-speed injection and multi-analyte processing which increases the productivity and throughput of your LC-MS/MS system. The exceptional injection precision and minimal carryover enables flawless performance for LC-MS/MS analysis. The sample capacity of the base QS-Works Autosampler bundle can be increased by including an additional three-drawer Peltier stack to meet your lab requirements.

## Hardware Specifications

Injection volume	Programmable from 0.1 $\mu$ L to 10000 $\mu$ L (depending upon loop and syringe size)
Injection Precision	Full Loop Injection: RSD $\leq$ 0.1 % for 2 $\mu$ L & 20 $\mu$ L (caffeine in water) Partial Loop Injection: RSD $\leq$ 0.5 % for 2 $\mu$ L & 20 $\mu$ L (caffeine in water)
Injection Linearity	R2 $\geq$ 0.9999
Needle wash	Fully programmable volumes and speed for inside and outside needle washes
Wash Station	2 different wash solvents
Valve Switching Time	100 msec
Injection Time (minimum)	< 20 sec. (injection to injection including needle wash)
Loop Volume	2 $\mu$ L to 5 mL, 10 $\mu$ L standard
Syringe Volume	Standard: 100 $\mu$ L, Other syringe sizes available
Pressure Limit	Up to 20,000 psi
CarryOver	$\leq$ 0.004 % with LCMS-P Tool; 100 $\mu$ L
Sample capacity	One three-drawer Peltier Stack supports up to: - 6 x 54 vial tray (standard 2 mL vials): 324 - 6 x 70 vial tray (1 mL vials): 420 - 6 x 15 vial tray (10 mL vials): 90 - 6 MT/DW plates <i>*An additional three-drawer Peltier Stack can be purchased to increase the sample capacity by two times.</i>

Wetted Parts in Flow Path PEEK, PFA, FEP, FFPM

Peltier Temperature Range 4 to 40°C

## Operating Specifications

Working Temperature	5 to 40°C (41 to 104°F)
Storage Temperature	-25 to +65°C (-13 to 149°F)
Humidity	Max. 80% RH, non-condensing
Work Environment	Indoor use
Altitude	Up to 3000 m above sea level
Sound Pressure Level	<70 dBA

## Physical Specifications

Power Requirements	100 - 240 VAC +/- 10%, 50 - 60 Hz
Dimensions	Width (X-Axis) x Depth (Y-Axis) x Height (Z-Axis) 53.4 x 50.3 x 54.7 cm (21.0 x 19.8 x 21.5 in.)
Weight	12.24 kg (26.98 lbs) without Peltier Stack 22.27 kg (49.10 lbs) with one three-drawer Peltier Stack

For more information  
visit [www.perkinelmer.com](http://www.perkinelmer.com)

PerkinElmer, Inc.  
940 Winter Street  
Waltham, MA 02451 USA  
P: (800) 762-4000 or  
(+1) 203-925-4602  
[www.perkinelmer.com](http://www.perkinelmer.com)

  
**PerkinElmer**<sup>®</sup>  
For the Better

For a complete listing of our global offices, visit [www.perkinelmer.com/ContactUs](http://www.perkinelmer.com/ContactUs)

Copyright ©2020, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.  
FOOD\_SPC\_20201113\_1.1 PKI