

## iQue® QMax - Extend Sheath and Waste Fluidic Run-Time

QMax Workstation takes fluid management off your plate.

- Automatically maintains sheath and waste levels in iQue Forecyt® software
- Nozzle emptying eliminates heavy waste container lifting
- Mobile trolley design for easy refill and waste removal
- Fully integrated with the automation interface (API)

QMax Workstation manages sheath and waste levels, so you don't have to.

- Extend hands-off runtime on the iQue® HTS platform by up to 5 fold \*
- Keep experiments running continuously with fluids that can be replaced while the system is running
- Increase productivity with additional plates or longer sampling times

\*Your runtime may vary.



# What's New in iQue® QMax?

- Extended sheath and waste runtime – up to 48 hours
- Quality-of-life improvements and safer maneuverability
- Enhanced automation with improved serial/TCP connectivity and software upgrades



*The new trolley is fantastic and we use the nozzle for easy waste removal everytime.*

**-iQue® User**



iQue® QMax Refill Module

iQue® QMax Workstation

	iQue® QMax Refill Module	iQue® QMax Workstation
Order No.	na	BA-97146
Total Sheath Capacity	20L	20L
Waste Sheath Capacity	20L	20L
iQue® Screener, Plus, 3 Compatible	Yes	Yes
iQue® 5 Compatible	Yes	Yes
Direct Connection to iQue Forecyt®	No	Yes
Mobile - Trolley	No	Yes
Easy Waste Clearance with Nozzle	No	Yes
QMax Refill Module Sheath Carboy Filter	#91311	#91311
Sheath Additive Concentrate (100 mL)	#91310	#91310

For More Information, Please Contact Your Sales Representative