

Calculate the Real Cost of a Pre-Owned Liquid Handler.

The full cost of ownership can catch you off guard, especially if you're about to buy an older robot. Here's a handy worksheet and example to help estimate your total investment. Use this to get a clear picture of what you're signing up for.

Item	Cost	Comment
Liquid handler		Be sure to compare apples to apples, is the available configuration what you need?
Software update		Many software licenses are operating system dependent, and some are non-transferable.
Work-deck reconfiguration		Changing a used system into your needed solution may demand repurposing with new carriers, devices and even arms.
Service contract		Many used providers have no warranty or reduced warranty periods.
Hardware updates due to wear & tear		Wear and tear issue increases with instrument age, as does the risk of interrupted runs and unplanned down time.
Application support		Expert support is needed to correctly configure the application that you're running on the instrument.
Admin cost, incl. shipping and setup		The eye-catching prices from an auction house or used equipment broker is often just for the instrument.
Cost of a failed run		Count in the cost of reagents, precious sample, and your reputation when things go wrong.

Need some help with the numbers for your specific situation?

Want to know the service history of the instrument you're considering buying?

Scan the QR code to have an expert contact you to walk through your worksheet and give you the history of the instrument.

