

# THE PERFECT WAY TO KEEP YOUR COOL



## Chillers to Keep Your Lab Cool

The **DuraChill® recirculating chiller** provides cooling power for demanding applications and serves as an economical alternative to tap water cooling systems. Easy to use and maintain, it combines technological innovation with precise temperature control to deliver reliable heat removal for a wide variety of applications.

Its intuitive color touchscreen display offers a choice of five operating languages, and its home screen simultaneously displays set temperature, actual temperature, fluid pressure, fluid level, air filter status, and timer.

At the same time, continuous fluid-level monitoring protects the pump and prevents freezing, and a 1 HP turbine pump provides reliable flow. What's more, its Whispercool® noise reduction system reduces operating noise levels to almost imperceptible levels while maintaining optimal cooling performance, even in high ambient-temperature operating conditions.



**DuraChill Chiller**

### **Additional features and benefits include:**

- Highly optimized design reduces refrigerant gas needed and allows for migration to natural refrigerants
- Intelligent self-diagnostic system compares the chiller's functionality to its original factory operating state to minimize performance issues
- Front-fill reservoir access
- Externally accessible adjustable pressure regulator valve
- Self-changing air filter system eliminates preventative maintenance by automatically changing the filter media at user-definable intervals (optional)
- UV Light Biological Growth Inhibitor System eliminates the need for ecosystem-damaging algacides (optional)



### DuraChill Dynamic Filter and UV Light Biological Inhibitor Option Part Number N0772052

- DuraChill is pre-programmed to change its filter once a month for a two-year period and can be adjusted to suit your own operating environment. This will reduce unnecessary service calls on your new chiller.
- As part of our commitment to the environment, we offer a UV light biological growth inhibitor that will continuously control biological growth in the fluid path without the need to add chemical growth inhibitors and prevent algicides from being released into the world's waterways.

**Dynamic Filter and UV Light Biological Inhibitor**

	DuraChill Chiller 240 V/50 Hz	DuraChill Chiller 230 V/60 Hz
<b>Part Number</b>	<b>N0772050</b>	<b>N0772051</b>
Power Requirements	240 V, 50 Hz 13.5 A	230 V, 60 Hz, 13.5 A
Cooling Capacity @20 °C	2650 Watts (9040 Btu/hr)	2900 Watts (9895 Btu/hr)
Cooling Capacity @10 °C	1900 Watts (6480 Btu/hr)	1930 Watts (6585 Btu/hr)
Cooling Capacity @0 °C	1000 Watts (3410 Btu/hr)	1000 Watts (3410 Btu/hr)
Pump Pressure Range	20 to 75 psi/138 to 621 kPa	20 to 90 psi/138 to 689 kPa
Maximum Pump Flow Rate	2.9 gpm/ 11 L/min	3.5 gpm/ 13.2 L/min
Fluid Temperature Range	-10 to 70 °C (14 to 158 °F)	
Temperature Stability	+/-0.1 °C (+/-1.8 °F)	
Display Temperature Units	°C or °F	
Display Pressure Units	psi or kPa	
Pump Inlet and Outlet	½ inch female NPT	
Connectivity	RS-232, USB Serial Port Emulation, USB TMC, Dry Contact On/Off/Status	
Selectable Operating Languages	English, Spanish, French, German, Mandarin	
Ambient Air Temperature Tracking	Yes	
Self Changing Air Filter	Optional	
UV Biological Growth Inhibitor	Optional	

**Shop Online Today!**  
<https://www.perkinelmer.com/buy>

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)



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.