

Site Requirements for the ICELL8® Multi-Sample NanoDispenser Plus

D14-000020-0001

Revision A

Date Effective: 07/14/2022

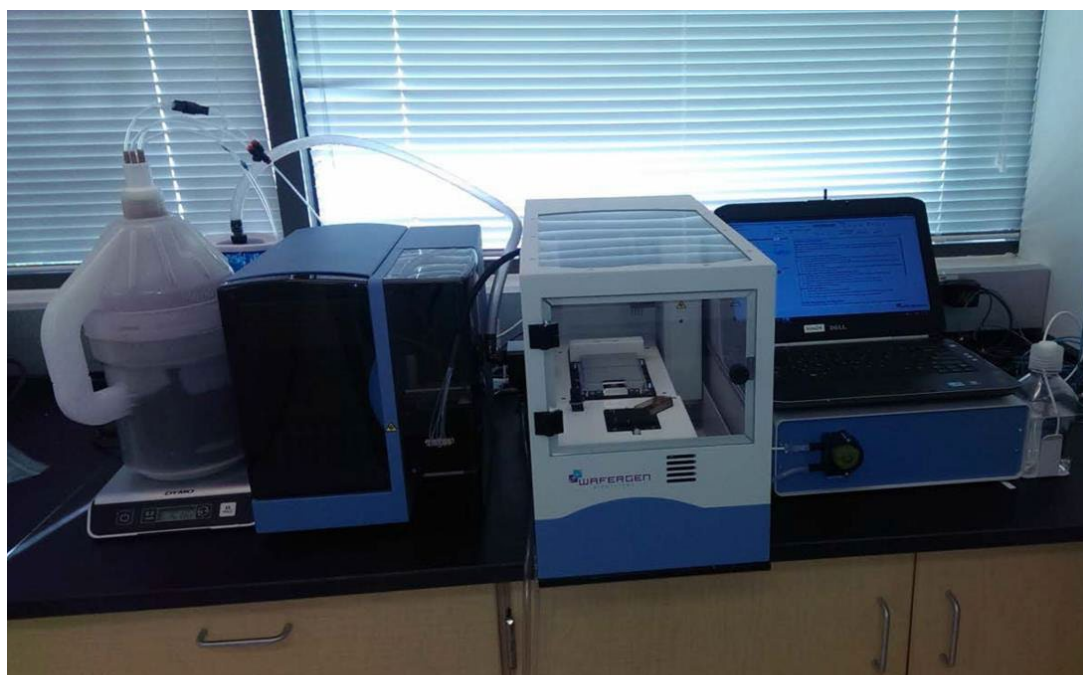
Parameter	Requirement	*Customer Initials	*Takara Bio Technician Initials
Bench space including clearance (capable of supporting approx. 50 kg, including CPU)	<ol style="list-style-type: none"> 1. Dispenser, Pump Box, CPU, Pressure reservoir, transformer (if required): 135 cm Wide x 75 cm Deep x 46 cm High 2. Humidifier 30.5 cm H x 20.5 cm W x 17.8 cm L 		
Floor space	<p><u>Sufficient to accommodate</u></p> <ol style="list-style-type: none"> 1. Helium (preferred) or argon source, 25 cm diameter bottle x approx. 152 cm tall 2. Waste Bottle 25 cm diameter x 46 cm tall 		
Power	<ol style="list-style-type: none"> 1. Power (either a OR b OR c) <ol style="list-style-type: none"> a. 120 VAC / 60 Hz mains One or more 20A circuits to power and a minimum of 4 ea. NEMA 5-15 receptacles for the following loads: <ol style="list-style-type: none"> i. Dispenser ii. Pump Box iii. Laptop computer (The entire system may be powered from a single 120V- 20A circuit). b. 230 VAC AC mains One 10A circuit with Schuko receptacles to power A 230:115/120 step-down transformer with provision for 4 ea. NEMA 5-15 receptacles (this transformer is available from Takara Bio) but requires a power strip with a minimum of three receptacles. The receptacles power the Dispenser, Pump Box, and Laptop Computer. c. 100 VAC / 50, 100V /60 Hz mains Two 15A circuits to power and a minimum of 4 ea. NEMA 5-15 receptacles for the following loads: <ol style="list-style-type: none"> i. Dispenser ii. Pump Box iii. Laptop computer 		

*By initialing, you acknowledge that you are verifying the site requirements set forth by this document.

ICELL8® Site Requirements: Multi-Sample NanoDispenser Plus

Parameter	Requirement	*Customer Initials	*Takara Bio Technician Initials
Helium (or Argon)	<ol style="list-style-type: none"> 99.9% purity or greater. Capacity Approximately 223 standard cubic feet (reported at 15.6°C and 1 atmosphere (1.01325 Bar). This capacity is sufficient for six months or more of typical usage Pressure: 2265 psi (150 Bar), with regulation from 0 to 200 psig (0-15 Bar). Typical usage pressure 30 psig (2 Bar) Fittings to accommodate 1/8" (3.2 mm) outer diameter flexible urethane tubing fittings (push-to-connect fittings). Acceptable thread forms are 1/8" NPT (female) or M5 straight thread (female). Consult Takara Bio for other thread forms. 		
Reagents	<ol style="list-style-type: none"> Filtered water: Milli-Q or Elga system Reagent grade 100% Isopropyl Alcohol for priming system 0.2% bleach solution (diluted from commercial bleach solution with a date of manufacture of less than a year in filtered water) 		
Environmental conditions	<ol style="list-style-type: none"> 15-30 °C Ambient 25-70% Relative Humidity, non-condensing <2000 meters altitude Pollution degree 2 or less 		
Information Technology Requirements	Customer's IT representative has read and understands Takara Bio document D06-000008 for IT requirements of SmartChip® Systems		

*By initialing, you acknowledge that you are verifying the site requirements set forth by this document.



Contact Us	
Customer Service/Ordering	Technical Support
tel: 800.662.2566 (toll-free)	tel: 800.662.2566 (toll-free)
fax: 800.424.1350 (toll-free)	fax: 800.424.1350 (toll-free)
web: takarabio.com/service	web: takarabio.com/support
e-mail: ordersUS@takarabio.com	e-mail: technical_support@takarabio.com

Notice to Purchaser

Our products are to be used for **Research Use Only**. They may not be used for any other purpose, including, but not limited to, use in humans, therapeutic or diagnostic use, or commercial use of any kind. Our products may not be transferred to third parties, resold, modified for resale, or used to manufacture commercial products or to provide a service to third parties without our prior written approval.

Your use of this product is also subject to compliance with any applicable licensing requirements described on the product's web page at takarabio.com. It is your responsibility to review, understand and adhere to any restrictions imposed by such statements

© 2017 Takara Bio Inc. All Rights Reserved.

All trademarks are the property of Takara Bio Inc. or its affiliate(s) in the U.S. and/or other countries or their respective owners. Certain trademarks may not be registered in all jurisdictions. Additional product, intellectual property, and restricted use information is available at takarabio.com.

This document has been reviewed and approved by the Quality Department.