

Specification: Site Requirements for Imaging System

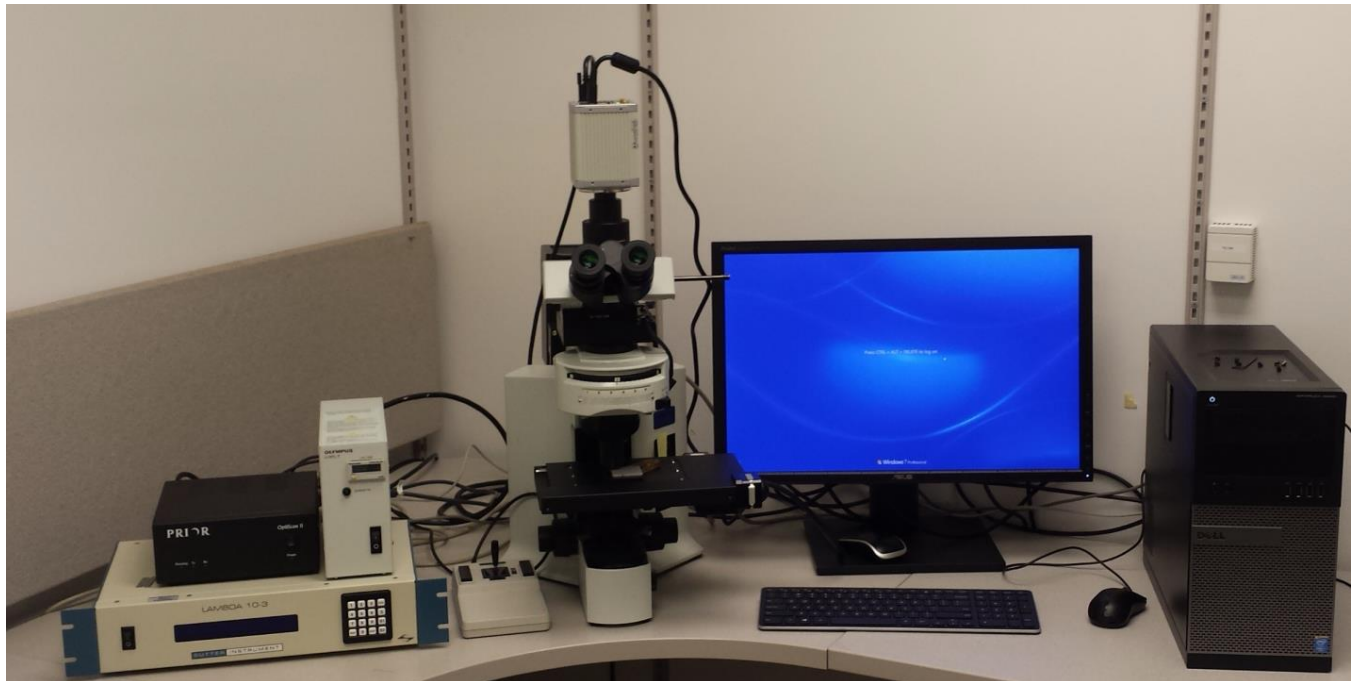
D14-000029-0001

Revision A

Date Effective: 05/18/2015

Parameter	Requirement	*Customer Initials	*Takara Bio Technician Initials
Bench Space (includes clearance distances).	Width: 64" Depth: 26" Height: 34"		
AC Power	100–240 VAC, 2.0/1.0 A (47–63 Hz), power outlet within 2 m Main supply voltage with ≤10% variance from nominal		
Environmental Conditions	15-30 °C Ambient 10-70% Relative Humidity, non-condensing Pollution degree 2 or less		
Optical	No direct sun light.		

*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	web: takarabio.com
e-mail: ordersUS@takarabio.com	e-mail: techUS@takarabio.com

Notice to Purchaser

Our products are to be used for research purposes only. They may not be used for any other purpose, including, but not limited to, use in drugs, *in vitro* diagnostic purposes, therapeutics, or in humans. 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 prior written approval of Takara Bio USA, Inc.

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.