

GeneTitan™ Multi-Channel Instrument

Catalog Numbers 00-0372 and 00-0373

Pub. No. MAN0025663 Rev. A.0

Customer Concierge contact information

Note: If your site or region is not covered by customer concierge service, contact your local Thermo Fisher Scientific service and support representatives for site preparation and instrument installation.

We are here to help. Send your questions to:

Location	Concierge for your region
North America	AMER.Concierge@thermofisher.com
Europe, Middle East, and Africa	EMEA.Concierge@thermofisher.com
Asia-Pacific and Japan	APJ.Concierge@thermofisher.com

Note: Complete the *GeneTitan™ Multi-Channel Instrument Site Preparation Checklist*, Pub. No. MAN0025663, then click **Submit** at the bottom of the checklist to send the form to your Concierge team. If your region is not supported by one of the teams listed above, email the completed checklist to your field application scientist.

Site preparation

IMPORTANT! To facilitate a successful and timely installation, ensure that this site preparation checklist is complete with all of the requirements checked. When complete, use the Submit button at the end of this document to send the information to the Thermo Fisher Scientific Customer Concierge team for your region.

Requirement	For details, see specific sections of the site preparation guide ^[1]
General	
<input type="checkbox"/> Customer responsibilities have been reviewed.	"Customer responsibilities"
<input type="checkbox"/> Personnel have been assigned tasks and responsibilities.	
The installation site is identified and meets the following requirements:	
<input type="checkbox"/> Space and clearance	"Configured system dimensions and connections"
<input type="checkbox"/> Pneumatic requirements	"Pneumatic requirements"
<input type="checkbox"/> Laboratory workbench	"Laboratory workbench requirements"
<input type="checkbox"/> Clean dry air	"Clean dry air (CDA)"
<input type="checkbox"/> Environmental	"Exhaust and venting requirements"
<input type="checkbox"/> Electrical	"Electrical outlets"
<input type="checkbox"/> Network	"Network requirements"
<input type="checkbox"/> Safety	"Safety requirements"

Requirement	For details, see specific sections of the site preparation guide ^[1]
Workstation	
The workstation will be connected to a network. <input type="radio"/> Yes <input type="radio"/> No (recommended) The pre-installed antivirus software on the workstation will be removed and replaced. <input type="radio"/> Yes <input type="radio"/> No If "Yes", provide the name of the replacement software. If "No", enter "N/A". <input type="text"/>	"Network requirements"
Receiving	
Do you have a loading dock? <input type="radio"/> Yes <input type="radio"/> No If "No", what special handling will you require? If "Yes", enter "N/A". <input type="text"/>	"Site requirements"
<input type="checkbox"/> The installation site is cleared and ready for the installation.	"Site requirements"
<input type="checkbox"/> All required customer-supplied materials for installation, qualification, and operation are available.	"Required materials not supplied"

^[1] GeneTitan™ Multi-Channel Instrument Site Preparation Guide, Pub. No. MAN0025571.

Customer information

Company name

Region North America
 Europe, Middle East, and Africa
 Asia Pacific and Japan
 Other

Installation site address

Facilities

Name

Phone

Email

IT personnel

Name

Phone

Email

Scientist or primary end user

Name

Phone

Email

Acknowledgement

I verify that the above information is correct and that the site will be prepared before the arrival of the Thermo Fisher Scientific service representative.

I confirm that the necessary IT, facilities personnel, and scientists (end users) will be available to work with Thermo Fisher Scientific staff during the week of installation.

Name:

Email:

Date:



Affymetrix Inc. | 3450 Central Expressway | Santa Clara, California 95051 USA

For descriptions of symbols on product labels or product documents, go to [thermofisher.com/symbols-definition](https://www.thermofisher.com/symbols-definition).

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Revision history: Pub. No. MAN0025663

Revision	Date	Description
A.0	05 Jan 2022	New document for GeneTitan™ MC Instrument site preparation.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2022 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.