

revvity

# chemagic 360 Instrument Order Guide



For research use only. Not for use in diagnostic procedures.

## Table of contents

**Introduction** ..... 2

**chemagic 360 Instrument** ..... 3

**Required Accessories**..... 4

**Optional Items and Services** ..... 5

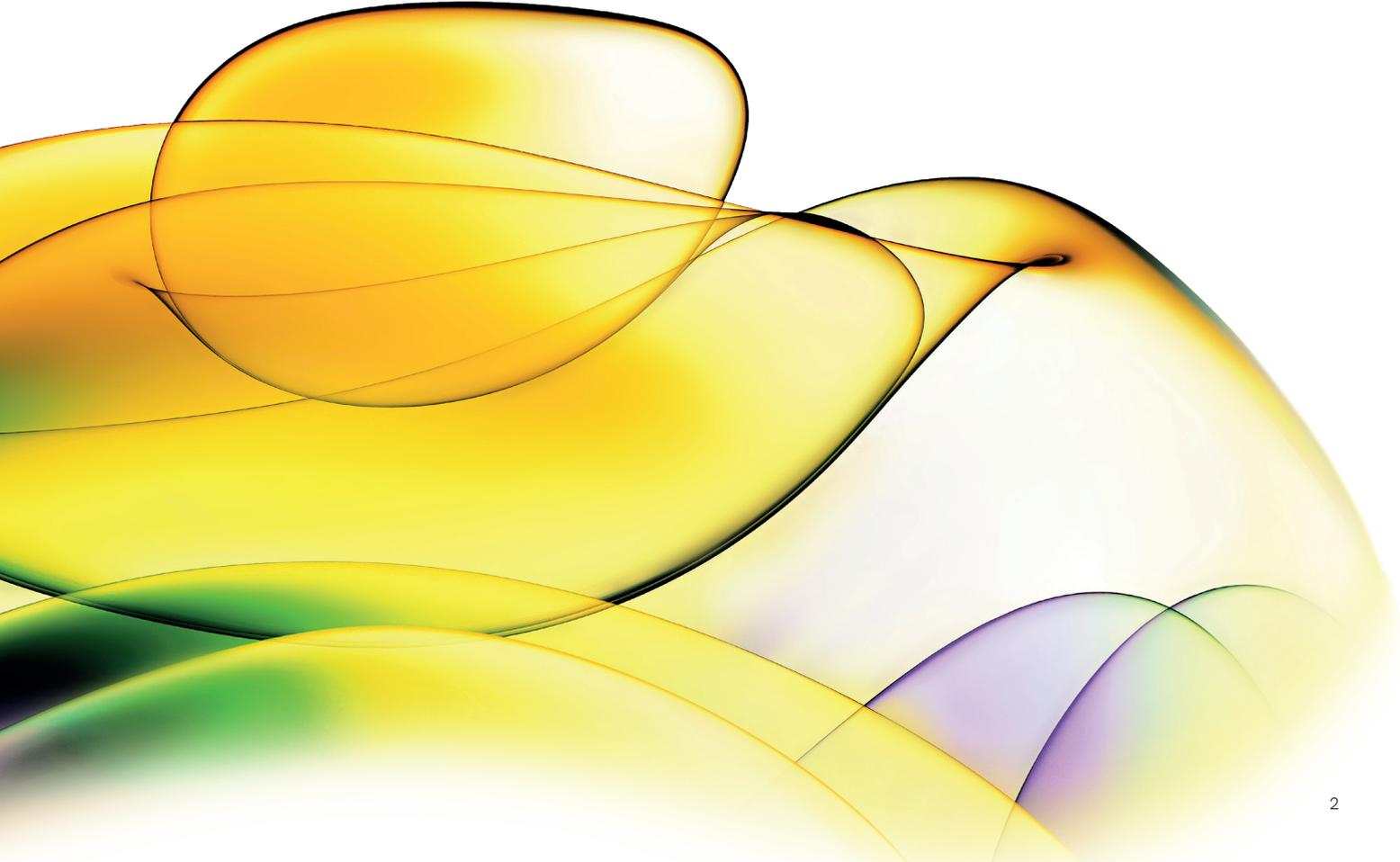
**Service Plans and Warranties** ..... 9

## Introduction

The chemagic™ 360 instrument is a trusted laboratory workhorse for quality automated nucleic acid purification from diverse sample materials. With easily interchangeable rod head sets, the system provides high flexibility, processing 12, 24, or 96 samples per run from a wide range of sample volumes. The instrument achieves high yields, purity and, integrity of nucleic acids based on reliable chemagic separation technology and chemagic kits that incorporate patented M-PVA Magnetic Beads.

This order guide summarizes the parts and services that are included or can be ordered in addition to your chemagic 360 instrument. Explore the range of options available to facilitate your workflow and learn about the services we can provide to promote smooth and effortless installation and use.

For any enquiries, please contact [support.chemagen@revvity.com](mailto:support.chemagen@revvity.com).





## chemagic 360 Instrument

The chemagic 360 instrument is sold as a stand-alone instrument. If you require the instrument to be integrated within an automated robotic workstation, please contact us for details.

Product number	Product name
2024-0020	<p><b>chemagic 360 Instrument</b></p> <p>The chemagic 360 instrument (RUO) is delivered with the following items:</p> <ul style="list-style-type: none"> <li>▪ Integrated PC which has a Windows Operating System (v10) in English and chemagic software in English</li> <li>▪ Keyboard</li> <li>▪ Monitor with touch screen display</li> <li>▪ EU standard cable (other cables are available on request)</li> <li>▪ Handheld barcode scanner</li> </ul>

For research use only. Not for use in diagnostic procedures.

## Required Accessories

The following items and services are not included with the instrument and must be ordered in addition.

### Installation Services

Product number	Service name and description
N0239033	<p><b>chemagic 360 IQ/OQ + Installation</b></p> <p>This is a service provided by trained personnel to ensure proper installation and operation qualification of the chemagic 360 instrument. This covers pre-installation checks, installation of equipment and software, and operational tests to certify standards of instrument functionality and communication are met.</p>

### chemagic Rod Head Sets

The chemagic Rod Head Sets include the chemagic Rod Heads with magnetizable steel rods that rotate using stirring motors, together with instrument accessories needed to support the 12-, 24-, or 96- throughput workflow. To cater to flexible sample input volumes and throughputs, chemagic Rod Heads are easily interchangeable. Please order at least one chemagic rod head set to be used together with the chemagic 360 instrument.



Product number	Product name																								
CMG-371	<p><b>chemagic 360 12 Rod Head Set</b></p> <p>The following components are included but can be ordered separately if required:</p>																								
	<table border="1"> <thead> <tr> <th>Part number</th> <th>Part name</th> </tr> </thead> <tbody> <tr> <td>CMG-534</td> <td>chemagic 12 Rod Head</td> </tr> <tr> <td>61010786</td> <td>Calibration Plate 12</td> </tr> <tr> <td>61010717</td> <td>Drop Cover Plate 360/12</td> </tr> <tr> <td>CMG-374</td> <td>Installation Kit 360/12</td> </tr> <tr> <td>CMG-375</td> <td>Manifold Unit 360/12</td> </tr> <tr> <td>CMG-510</td> <td>chemagic Rack Set 12</td> </tr> <tr> <td>CMG-308</td> <td>chemagic Stand 12</td> </tr> <tr> <td>CMG-594</td> <td>chemagic Lid-Set 6</td> </tr> <tr> <td>CMG-600</td> <td>chemagic Canister Stand for chemagic Dispenser</td> </tr> <tr> <td>CMG-13005116</td> <td>Waste Canister with quick coupling</td> </tr> <tr> <td>CMG-572</td> <td>chemagic Cleaning Bottle</td> </tr> </tbody> </table>	Part number	Part name	CMG-534	chemagic 12 Rod Head	61010786	Calibration Plate 12	61010717	Drop Cover Plate 360/12	CMG-374	Installation Kit 360/12	CMG-375	Manifold Unit 360/12	CMG-510	chemagic Rack Set 12	CMG-308	chemagic Stand 12	CMG-594	chemagic Lid-Set 6	CMG-600	chemagic Canister Stand for chemagic Dispenser	CMG-13005116	Waste Canister with quick coupling	CMG-572	chemagic Cleaning Bottle
	Part number	Part name																							
	CMG-534	chemagic 12 Rod Head																							
	61010786	Calibration Plate 12																							
	61010717	Drop Cover Plate 360/12																							
	CMG-374	Installation Kit 360/12																							
	CMG-375	Manifold Unit 360/12																							
	CMG-510	chemagic Rack Set 12																							
	CMG-308	chemagic Stand 12																							
	CMG-594	chemagic Lid-Set 6																							
	CMG-600	chemagic Canister Stand for chemagic Dispenser																							
	CMG-13005116	Waste Canister with quick coupling																							
	CMG-572	chemagic Cleaning Bottle																							



Product number	Product name	
CMG-376	<b>chemagic 360 24 Rod Head Set</b>	
	The following components are included but can be ordered separately if required:	
	Part number	Part name
	CMG-535	chemagic 24 Rod Head
	61010785	Calibration Plate 24
	61010718	Drop Cover Plate 360/24
	CMG-379	Installation Kit 360/24
	CMG-369	Manifold Unit 360/96/24
	CMG-1783	chemagic Rack set 24
	CMG-308	chemagic Stand 12
	CMG-594	chemagic Lid-Set 6
CMG-600	chemagic Canister Stand for chemagic Dispenser	
CMG-13005116	Waste Canister with quick coupling	
CMG-572	chemagic Cleaning Bottle	



Product number	Product name	
CMG-370	<b>chemagic 360 96 Rod Head Set</b>	
	The following components are included but can be ordered separately if required:	
	Part number	Part name
	CMG-536	chemagic 96 Rod Head
	61010784	Calibration Plate 96
	61010718	Drop Cover Plate 360/24
	11320170	96 square well plate
	CMG-369	Manifold Unit 360/96/24
	CMG-501	chemagic Rack Set 360/96
	11105245	Disposable tip rack, 96 tips
	CMG-301	chemagic Stand 96
	CMG-594	chemagic Lid-Set 6
	CMG-600	chemagic Canister Stand for chemagic Dispenser
	61011358	Reagent Bottle Unit
CMG-13005116	Waste Canister with quick coupling	
CMG-572	chemagic Cleaning Bottle	

## Optional Items and Services

To facilitate your workflow, these items can be installed on your instrument and allow you to incorporate a range of features such as heated lysis, UV disinfection, multi-kit use and more. In addition, services are available for the on-site setup or support of custom applications.



### Product number

61014370

### Product name and description

#### Device Table chemagic 360

This is a table that can be used to hold the chemagic 360 instrument. It has dimensions of 90 x 82 x 77 cm (W x H x D) and comes with a storage cabinet and extendable surface for the placement of buffer bottles. Shipping dimensions and weight are 100 x 100 x 120 cm, 80 kg.



CMG-429

#### chemagic 360 UV Disinfection Option

This item allows for decontamination of any residual nucleic acids using a compact airflow exchange chamber with a built-in UV lamp, fan, dust filters and control unit. Safety to the user is ensured while active airflow provides efficient decontamination that can be controlled by protocols adapted to the user's need.



CMG-13005658

#### chemagic 360 Portrait Dispenser Kit

This item allows for the dispensing of buffers in portrait format with 8 nozzles instead of landscape format with 12 nozzles. This may facilitate the running of partial plates or to match dispensing formats of upstream or downstream liquid handlers.



CMG-13009700

#### chemagic Heating Rack

The chemagic Heating Rack is a device used to heat up 96-deep well plates. This can be installed in one position on the chemagic 360 to enable heated lysis or elution. Temperatures can be set up to 100 degrees Celsius.



CMG-13009703

#### chemagic Turntable Power Adapter

This device allows for electrical output to two positions on the chemagic 360 instrument turntable. This enables two accessories that require power (e.g., heating rack, sleeve clamp rack) to be installed simultaneously with an uninterrupted power supply.



Product number	Product name and description
CMG-424	<p><b>chemagic Switchbox</b></p> <p>The chemagic Switchbox enables two kits to be connected to the chemagic 360 instrument for easy switching between kits and respective extraction workflows. Automatic selection of the kit can be done with the protocol or by rod head recognition.</p>
CMG-13009704	<p><b>chemagic 360 8-Pump Landscape Dispenser Kit</b></p> <p>This item consists of 8 pumps (2 additional) and 2 manifolds with 3 inlets to allow for in process cleaning or prefilling of additional reagents. Dispensing takes place in landscape format with 12 nozzles.</p>
CMG-13009705	<p><b>chemagic 360 8-Pump Portrait Dispenser Kit</b></p> <p>This item consists of 8 pumps (2 additional) and 2 manifolds with 3 inlets to allow for in process cleaning or prefilling of additional reagents. Dispensing takes place in portrait format with 8 nozzles.</p>
CMG-13009706	<p><b>chemagic 360 Interior Light and Window</b></p> <p>This item consists of an LED light situated within the chemagic 360 that allows for clearer visualization of the manifold and dispensing system. A transparent side window is also included to enable observation from outside the instrument.</p>
CMG-13009707	<p><b>chemagic 360 Status Light</b></p> <p>This is a USB interface LED Tower light that connects via USB port to the computer and is pole mounted to extend 50 cm above the chemagic 360 instrument. It comprises of coloured LED lights to show instrument status.</p>
CMG-382	<p><b>chemagic Sleeve Clamp Rack 360/96</b></p> <p>The chemagic Sleeve Clamp Rack 360/96 is an electric clamp rack for chemagic Tip Rack 96 that serves to hold the tip rack in place while the rod head picks up the tip sleeves.</p>



Product number	Product name and description
 <p>CMG-13005670</p>	<p><b>DWP XL Frame Notched</b></p> <p>A framed adaptor which prevents the user from placing the deep well plate in the wrong orientation, ensuring well A1 will always be in the same position under the rod head. This is to be placed in a position on the round table to ensure proper sleeve removal for the 24 rod head.</p>
 <p>CMG-425</p>	<p><b>DWP Adapter Notched</b></p> <p>An adaptor which prevents the user from placing the deep well plate in the wrong orientation, ensuring well A1 will always be in the same position under the rod head. This is placed on one position of the round table for the 96 rod head.</p>
 <p>CMG-13005660</p>	<p><b>DWP XL Adapter Notched</b></p> <p>An adaptor which prevents the user from placing the deep well plate in the wrong orientation, ensuring well A1 will always be in the same position under the rod head. This is placed on one position of the round table for the 24 rod head.</p>
 <p>CMG-13005340</p>	<p><b>chemagic 13ml double-bottom rack H24</b></p> <p>This item is required for cfDNA extraction workflows on the chemagic 360 equipped with the chemagic 24 Rod Head set. Three quantities are needed when running the workflow and have to be ordered separately.</p>
 <p>CMG-425</p>	<p><b>Tip Tray 96 Adapter</b></p> <p>This item enables the use of hanging tip trays which hold the sleeves that cover the magnetic rods. This is the preferred sustainable alternative to having sleeves delivered in plastic tip boxes.</p>
 <p>CMG-1305653</p>	<p><b>Lid Opener</b></p> <p>This item is used for loosening caps of reagent bottles delivered in chemagic extraction kits.</p>

## Optional Services

Product number	Service name and description
CMG-13005179	<p><b>1 Day Application Support</b></p> <p>The includes one day of on-site applications support for the development of user applications on the chemagic 360 instrument. It may include standard instrument familiarization and/or training.</p>

## Service Plans and Warranties

Service contracts offer dependable instrument repair and maintenance to minimize downtime, extending the life of your instrument. Choose between the three tiers of service plans for your chemagic 360 instrument:

**Platinum** - For laboratories where instrument uptime is critical to performance. Includes a proactive alert system, for enabled instruments, that triggers corrective actions to help avoid equipment downtime and unnecessary expenses.

**Gold** - Enhance laboratory productivity by maintaining instrument efficiency; with a priority breakdown response combined with software and technical support.

**Silver** - For laboratories looking to optimize instrument performance, regular preventative maintenance (PM) is essential.

A one year warranty is included with your chemagic 360 instrument but extended warranties are also available upon purchase, please refer to your local Territory Account Manager for more details.

Product number	Product name
NAP2024-0020PPT1	Platinum Plan 1PM - chemagic 360
NAP2024-0020PAU1	Gold Plan 1PM - chemagic 360
NAP2024-0020PAG1	Silver Plan 1PM - chemagic 360

Product number	Product name
NAP2024-0020PPT2	Platinum Plan 2PM - chemagic 360
NAP2024-0020PAU2	Gold Plan 2PM - chemagic 360
NAP2024-0020PAG2	Silver Plan 2PM - chemagic 360

Service contracts may differ across territories, please refer to your local Territory Account Manager for more details.



[www.revivity.com](http://www.revivity.com)

revivity

Revvity, Inc.  
940 Winter Street  
Waltham, MA 02451 USA  
[www.revivity.com](http://www.revivity.com)

For a complete listing of our global offices, visit [www.revivity.com](http://www.revivity.com)  
Copyright ©2023, Revvity, Inc. All rights reserved.

AG012312\_07\_OG