

KingFisher™ Apex Dx Heads

KingFisher™ Apex Dx 96 PCR Head, KingFisher™ Apex Dx 96 Combi Head, KingFisher™ Apex Dx 96 Deep Well Head, and KingFisher™ Apex Dx 24 Combi Head

Catalog Numbers B24079910, B24079920, B24079930, B24079940

Pub. No. MAN0029197 Rev. A.0



WARNING! Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from [thermofisher.com/support](https://www.thermofisher.com/support).

Product description

The KingFisher™ Apex Dx Heads are used on the KingFisher™ Apex Dx Purification Instrument only. The instrument allows changing of the magnetic heads in order to use different plates for sample processing.

For information about changing the heads see, KingFisher™ Apex Dx Purification Instrument Instructions for Use (IVD Workflow) (Pub. No. MAN0026509)

Intended use

The KingFisher™ Apex Dx Heads are intended for *in vitro* diagnostics (IVD) use as accessories for the KingFisher™ Apex Dx Purification Instrument to automate the process of isolation and purification of nucleic acids, proteins, and cells for IVD procedures in combination with a dedicated purification kit. The head is intended for use by professional users, such as technicians and physicians trained in molecular biological techniques.

Magnetic heads



WARNING! This product contains very strong permanent magnets. People wearing a pacemaker or metallic prosthesis should not use this product. A pacemaker or prosthesis may be affected or damaged if it comes in close contact with a strong magnetic field.

IMPORTANT! Keep magnetic heads away from each other and other magnets at all times. Clashing of the magnets can cause serious damage to the magnets.

- Do not place the magnetic head on top of the instrument or on any metal surfaces.
- Do not use metal tools when handling magnetic heads.
- Do not place the magnetic head in close proximity to magnetic tapes, computer discs or other magnetic storage systems, such as credit cards, as damage can occur due to the strong magnetic field of the magnetic heads.
- Do not place the magnetic heads near a PC display, as this may cause damage to the display.
- When not in use, keep the magnetic heads in their respective storage boxes.

Contents and storage















The KingFisher™ Apex Dx Purification Instrument is to be used together with one of four different magnetic heads detailed in the following table.

Contents ^[1]	Catalog No.
KingFisher™ Apex Dx 96 PCR Head	B24079910
KingFisher™ Apex Dx 96 Combi Head	B24079920
KingFisher™ Apex Dx 96 Deep Well Head	B24079930
KingFisher™ Apex Dx 24 Combi Head	B24079940

^[1] Store the magnetic heads in the provided box when they are not installed on the instrument.

Medical device symbols

The following table describes symbols that may be displayed on product labels. The symbols that are used on labels conform to standards *BS EN ISO 15223-1:2021* and *FDA 21 CFR 809.10 "Labeling for in vitro diagnostic products"*.

	IN VITRO DIAGNOSTIC MEDICAL DEVICE		MANUFACTURER
	CONSULT INSTRUCTIONS FOR USE		CATALOG NUMBER
	SERIAL NUMBER		AUTHORISED REPRESENTATIVE IN THE EUROPEAN COMMUNITY
	CAUTION		DATE OF MANUFACTURE
	CONTAINS		TEMPERATURE LIMIT
	UNIQUE DEVICE IDENTIFIER		INDICATES CONFORMITY WITH EUROPEAN UNION REQUIREMENTS
	INDICATES CONFORMITY WITH UNITED KINGDOM REQUIREMENTS		CAUTIOUS, STATIC MAGNETIC FIELD HAZARD

Serious incidents

Any serious incident that has occurred in relation to the device shall be reported to the manufacturer and the relevant regulatory authority in which the user and/or the patient is established.

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.



Life Technologies Holdings Pte Ltd | Block 33
| Marsiling Industrial Estate Road 3 | #07-06,
Singapore 739256



Life Technologies Europe B.V. Kwartsweg 2,
2665 NN Bleiswijk The Netherlands



Life Technologies Ltd 3 Fountain Drive
Inchinnan Business Park Paisley PA4
9RF Scotland, United Kingdom

Revision history: Pub. No. MAN0029197 A.0

Revision	Date	Description
A.0	9 May 2023	New document for KingFisher™ Apex Dx Heads.

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.

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.

©2023 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.