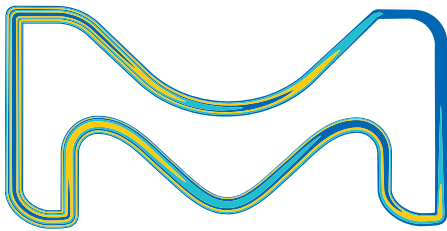


Application Pak Point-of-Use Polishers

Match water quality to your
every application



The life science business of
Merck KGaA, Darmstadt,
Germany operates as
MilliporeSigma in the
U.S. and Canada.

Milli-Q®
Lab Water Solutions

A range of Application Paks to fit your needs

Application Paks are specially designed point-of-use purifiers. When placed at the outlet of a POD water dispenser of a Milli-Q® IQ series water purification system, Application Pak point-of-use polishers ensure the water is free from contaminants that are critical to your applications.

Millipak® and sterile Millipak® Gold filters

A safe and efficient barrier against particulates and bacteria



Millipak® filters are designed for applications requiring the removal of particulates and bacteria. The filters are designed with a 0.22 µm PES (polyethersulfone) asymmetric membrane filter heat-sealed on a SAN (styrene acrylonitrile) housing. The membrane's conical pores allow high flow rate at low differential pressure; the membrane, housing material, and production processes have been selected to minimize organic and inorganic extractable release.

The Millipak® Gold 0.22 µm sterile filter has been gamma-irradiated for additional bacteria control.

Benefits

- Each Millipak® filter is individually tested and delivered with a Certificate of Quality
- Validated to deliver bacteria-free water (< 0.01 cfu/mL when installed and used in a laminar flow hood)
- Ease of maintenance: the filter is easily fitted and replaced

Biopak® Ultrafiltration Polisher

For pyrogen-, nuclease-, protease- and bacteria-free water



Designed for applications such as molecular biology, biochemistry, and cell culture, the Biopak® ultrafiltration (UF) cartridge removes macromolecules and larger biological structures such as bacteria. The device is made of polysulfone ultrafiltration hollow fibers. It is validated for the removal of pyrogens (endotoxins), RNases, DNases, proteases and bacteria.

The Biopak® ultrafiltration polisher provides a convenient and economical solution for scientists who need ultrapure water for cell culture, biochemistry or molecular biology applications.

Benefits

- Each Biopak® unit is delivered with a Certificate of Quality
- Validated to deliver pyrogen-free (< 0.001 EU/mL), RNase-free (< 1 pg/mL), DNase-free (< 5 pg/mL), protease-free (< 0.15 µg/mL) and bacteria-free water (< 0.01 cfu/mL)
- Eliminates the need for DEPC, therefore saving time, reducing costs, and removing risks
- Ease of maintenance: the polisher is easily fitted and replaced; does not require sanitization

Benefits to point-of-use purification with Application Pak polishers:

- **Versatility:** Adapt your water system to the various applications in your lab
- **Ease of use:** No maintenance required—simply replace the Application Pak when required by your water system; An ergonomic pack locker and protective bell provides 'snap-into-place' installation
- **Ensured data traceability:** e-Sure tags make it easy to automatically record Application Pak data and events throughout their lifetime

LC-Pak® Polisher

Water for trace and ultra-trace organic analyses



Designed for organic analyses by HPLC, UHPLC, LC-MS, and LC-MS/MS, the LC-Pak® Polisher is a device containing C18 reversed phase porous granular silica material. It is validated to achieve a water quality suitable for trace and ultra-trace analyses.

The LC-Pak® Polisher provides efficient and convenient access to freshly purified ultrapure water suitable for the most demanding trace and ultra-trace organic analyses.

Benefits

- Each LC-Pak® unit is delivered with a Certificate of Quality
- Provides water passing the following tests: HPLC gradient test at 210 and 254 nm; UV absorbance within specifications at 200 nm, 205 nm, 210 nm, and 254 nm; fluorescence as quinine < 1 ppb at 254 and 365 nm; Reserpine test: no peak > 10 ppb; residue after evaporation < 0.0001% w/w as described in ISO® 3696
- Delivers a minimum of 500 L of ultrapure water suitable for trace or ultra-trace organic analyses
- Ease of maintenance: the polisher is easily fitted and replaced

VOC-Pak® Polisher

Water for analysis of volatile organic compounds



Designed for the production of ultrapure water suitable for the analysis of volatile organic compounds (VOCs) in water by GC and GC-MS, the VOC-Pak® unit is a device made of activated carbon material specifically designed for efficient VOC removal.

The VOC-Pak® Polisher provides an efficient solution for scientists who need volatile organic compound-free ultrapure water for their laboratory work.

Benefits

- Each VOC-Pak® unit is delivered with a Certificate of Quality
- Provides water suitable for VOC analysis by gas chromatography
- Delivers a minimum of 300 L of VOC-free ultrapure water
- Ease of maintenance: the polisher is easily fitted and replaced

EDS-Pak® Polisher

Water for endocrine disruptor experiments

Designed for the production of endocrine disruptor-free ultrapure water, the EDS-Pak® polisher is a device made of a specific activated carbon material. The activated carbon used was selected for its specific and fast adsorption of endocrine disruptors. The EDS-Pak® polisher is validated to deliver water with a low level of endocrine disruptors, including phthalate esters, bisphenol A and nonylphenol.

The EDS-Pak® Polisher provides scientists with a convenient solution for their endocrine disruptor-free ultrapure water needs.

Benefits

- Each EDS-Pak® unit is delivered with a Certificate of Quality.
- Provides water suitable for endocrine disruptor analyses.
- Ease of maintenance: the polisher is easily fitted and replaced.



Ordering Information

Product	Cat. No.	Product	Cat. No.
Millipak® 0.22 µm Filter		LC-Pak® Polisher	
 Bacteria-free and particulate-free water	MPGP002A1	 Water for trace and ultra-trace organic analysis	LCPAK00A1
Millipak® Gold 0.22 µm Sterile Filter		VOC-Pak® Polisher	
 Bacteria-free and particulate-free water (gamma-irradiated)	MPGPG02A1	 Water for volatile organic compounds analysis	VOCPAK0A1
Biopak® Polisher		EDS-Pak® Polisher	
 Pyrogen-, nuclease-, protease- and bacteria-free water	CDUFBI0A1	 Water for endocrine disruptor experiments	EDSPAK0A1

To place an order or receive technical assistance

In the U.S. and Canada, call toll-free 1-800-645-5476
For other countries across Europe and the world,
please visit: [EMDMillipore.com/offices](https://www.emdmillipore.com/offices)
For Technical Service, please visit: [EMDMillipore.com/techservice](https://www.emdmillipore.com/techservice)

MilliporeSigma
400 Summit Drive
Burlington, MA 01803

[EMDMillipore.com](https://www.emdmillipore.com)

© 2019 Merck KGaA, Darmstadt, Germany and/or its affiliates. All Rights Reserved. MilliporeSigma, the vibrant M, Milli-Q, Millipak, EDS-Pak, Biopak, LC-Pak and VOC-Pak are trademarks of Merck KGaA, Darmstadt, Germany and/or its affiliates. All other trademarks are the property of their respective owners. Detailed information on trademarks is available via publicly accessible resources.

