



Simplifying your manual platform DELFIA™ workflow

DELFIA™ Trio modernizes current manual DELFIA workflows with intelligent instrument 'quality control' checks all supported by an intuitive touch screen designed to support ease of use in busy laboratory workflows.

The 1296-0070 DELFIA Trio is a 3-in-1 instrument developed for sample processing of DELFIA manual kits. It combines disk removing, washing, and dispensing into one instrument offering a streamlined workflow for manual DELFIA users.

DELFIA Trio has been designed to improve your workflows by:

Saving Your Precious Time

- We've designed DELFIA Trio to support your busy lab management practice. Providing you more time to focus on important day to day priorities.

'Load & Go' Simplification Runs

- Optimising your processing steps can now be found in ONE Instrument!
- Disk Removing - Washing - Dispensing, all combined and self-contained

Building Stronger Lab Management Practice

- Process Quality Checks in built
- Easy Read Intuitive Touchscreen
- 14 pre-installed DELFIA protocols



- ✔ Saves Your Precious Time
- ✔ 'Load & Go' Simplification Runs
- ✔ Stronger Lab Management Practice



Simple intuitive user interface

| Physical data | | |
|--------------------------------|-------------------|--|
| Dimensions | Width | 538 mm (instrument), 1102 mm (with bottles) |
| | Length | 590 mm |
| | Height | 336 mm (instrument), 527 mm (with bottles) |
| | Weight | Instrument 37 kg, ~ 50 kg (with bottles and bottle holders) |
| Operating conditions | Temperature | +15°C to +30°C |
| | Relative humidity | 10 % to 80 % relative humidity |
| Electrical requirements | Voltage/Frequency | 100 - 240 VAC / 50/60 Hz |
| | Consumption | 160 W |
| | Fuses | 2 x 3.15 A (24 V; 5,25 A) |
| | Power cord | External universal PSU, cord according to region |
| Hardware specifications | | |
| Manifold | | 12 channels |
| Volumes of bottles | Waste | 5 L |
| | Waste | 2 L |
| | Wash | 2 L |
| | Rinse | 2 L |
| User interface | | 7" touchscreen |
| Other features | | Bottle holder with a shield for enhancement solution |
| Software specifications | | |
| Protocols | | 14 ready-made factory protocols for manual DELFIA assays, possibility for custom protocols |
| Number of strips | | 2, 4, 6 or 8 |
| Disk removal | | Yes/no |
| No. of wash cycles | | 0 to 20 |
| Dispensing | | Yes/no (200 µl) |
| Plates | | Nunc 8x12 or Thermo 8x12 Immunostrips |
| Other features | | Pressure analysis based monitoring of protocol completion status |
| | | Maintenance wizard |
| | | Possibility to inactivate modules |
| Technical features | | |
| Vacuum data | | Flow rate in normal atmosphere 14.5 l/min, final vacuum 72 % |
| Switch positions | | Power switch on back panel, power button on front panel |

Products may not be licensed in accordance with the laws in all countries, such as the United States and Canada. Please check with your local representative for availability. Please note that product labeling (such as kit insert, product label, and kit box) may be different compared to the company branding. Please contact your local representative for further details.