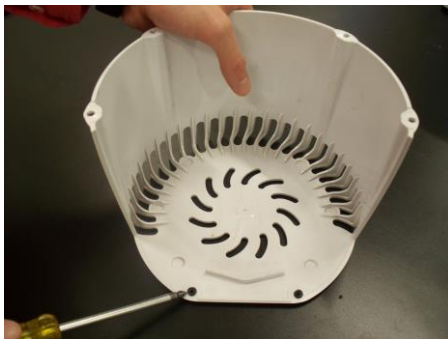
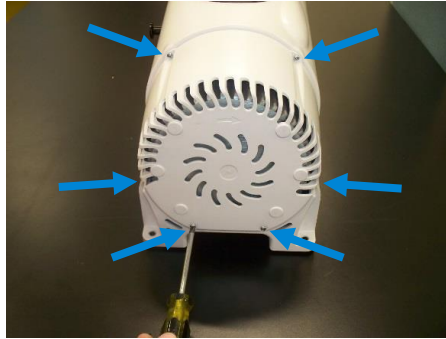


IDP7/IDP10 Rear Cowling Replacement Procedure



Equipment List:

- #2 Phillips screw driver

Disassembly:

WARNING



Turn off pump and unplug AC power cord. As some surfaces on the interior of the pump can become uncomfortably warm to the touch during operation, Agilent recommends the pump be allowed to cool to room temperature before beginning the operation.

- Remove (6) 4 mm screws from rear cowling.
- Discard (2) rear cowling rubber washers.

Assembly:

- Insert (2) 4 mm screws through the bottom (2) holes.
- Thread on rear cowling rubber washers onto the (2) screws, as shown.
- Position rear cowling to align bottom two screws with washers in to the base cowling.
- Thread the bottom (2) screws into base cowling, but not fully tightened.
- Thread the remaining (4) screws into the base cowling, but not fully tightened.

CAUTION



Cautions appear before instructions, which if not followed, could cause damage to the equipment or data loss.

WARNING



Warnings appear for a particular procedure or practice which, if not followed correctly, could lead to serious injury or death.

CAUTION



Do not overtighten screws.

- **Secure all (6) screws.**

United States
Agilent Technologies
121 Hartwell Avenue
Lexington, MA 02421 - USA
Tel. +1 781 861 7200
Fax. +1 781 860 5437
Toll-Free: +1 800 882 7426
vpl-customer-care@agilent.com

Belgium
Agilent Technologies Belgium S.A./N.V.
Customer Contact Center
Pegasus Park
De Kleetlaan 12A bus 12
B-1831 Diegem
Tel. +32 2 404 92 22
Fax. +32 2 626 46 30
customer-care_belgium@agilent.com

Brazil
Agilent Technologies Brasil
Avenida Marcos Penteado de Ulhoa
Rodrigues, 939 - 6º andar
Castelo Branco Office Park
Torre Jacarandá - Tamboré
Barueri, Sao Paulo CEP: 06460-040
Toll-Free: +0800 728 1405

China
Agilent Technologies (China) Co. Ltd
No.3, Wang Jing Bei Lu, Chao Yang District
Beijing, 100102, China
Tel. +86 (0)10 64397888
Fax. +86 (0)10 64392765
Toll free: 800 820 6778 (Phone)
Toll free: 400 820 6778 (Cell Phone)
vacuum.cnmarketing@agilent.com
vpc-customerservice@agilent.com

France
Agilent Technologies
Parc Technopolis - Z.A. de Courtaboeuf
3, avenue du Canada - CS 90263
91978 Les Ulis cedex, France
Tel. +33 (0) 1 64 53 61 15
Fax. +33 (0) 1 64 53 50 01
vpf.sales@agilent.com

Germany and Austria
Agilent Technologies
Sales & Services GmbH & Co. KG
Lyoner Str. 20
60 528 Frankfurt am Main
GERMANY
Ph. +49 69 6773 43 2230
Fax. +49 69 6773 43 2250

India (Sales)
Agilent Technologies India Pvt. Ltd.
Unit Nos 110-116, & Part of 101 & 109
First Floor, Splendor Forum,
Plot No. -3, District Centre, Jasola
New Delhi-110025
Tel. +91 11 4623 7100
Fax. +91 4623 7105

India (Service)
Agilent Technologies India Pvt. Ltd.
C-Block, RMZ Centennial Plot Number -8A,
8B, 8C, 8D,
Doddanakundi Industrial Area, ITPL Road,
Mahadevapura Post, Bangalore-560048
Tel. +91 80 40614000
Fax. +91 80 40148991

Italy
Agilent Technologies Italia S.p.A.
Via F.lli Varian, 54
10040 Leini, (Torino) - Italy
Tel. +39 011 9979 111
Fax. +39 011 9979 350
Toll free: 00800 234 234 00
vpt-customerservice@agilent.com

Japan
Agilent Technologies Japan, Ltd.
9-1 Takakura-cho Hachioji-city,
Tokyo, Japan
Tel. +81-3-5232-1253
Fax. +81-120-565-154
Toll-Free: +81-120-477-111
jp-vvt-sales.pdf-ext@agilent.com

Korea
Agilent Technologies Korea, Ltd.
Ilshin Building 4F
Yongsan-gu Hannam-daero
Seoul Korea 04418
Tel. +82 (0)2 2194 9449
Fax. +82 (0)2 2194 9853
Toll free: 080 222 2452
vpk-customerservice@agilent.com

Mexico
Agilent Technologies
Concepcion Beistegui No. 109
Col Del Valle
C.P. 03100 - Mexico, D.F.
Tel. +52 5 523 9465
Fax. +52 5 523 9472

Netherlands
Agilent Technologies Netherlands B.V.
Customer Contact Center
Laan van Langerhuize 1, toren A-8
1186 DS Amstelveen
Tel. +31 020 547 2600
Fax. +31 020 654 5748
customer-care_netherlands@agilent.com

Singapore
Agilent Technologies Singapore Pte. Ltd,
1 Yishun Avenue 7,
Singapore 768923
Tel. (65) 6215 8045
Fax. (65) 6754 0574
Toll-Free: 1 800 2762622
vps-customerservice@agilent.com

Southeast Asia
Agilent Technologies Sales Sdn Bhd
Unit 201, Level 2 uptown 2,
2 Jalan SS21/37, Damansara Uptown
47400 Petaling Jaya, Selangor, Malaysia
Tel. +603 7712 6181
Fax. +603 7727 1239
Toll-Free: +1 800 880 805
vps-customerservice@agilent.com

Taiwan
Agilent Technologies Taiwan Limited
No. 20 Gao-shuang Road, Ping-zhen Dis
Tao-Yuan City
32450 Taiwan, R.O.C.
Tel. +8863 3 4959004
Toll-Free: 0800 018 768
vpw-customerservice@agilent.com

UK and Ireland
Agilent Technologies LDA UK Limited
Lakeside Cheadle Royal Business Park
Cheadle, Cheshire SK8 3GR,
United Kingdom
Tel. +44 01865291570
Fax. +44 01865291571
Toll-Free: 00 800 234 234 00
Toll-Free fax.: 00 800 345 345 00
vpt-customer-care@agilent.com

Other Countries
Agilent Technologies Italia S.p.A.
Via F.lli Varian, 54
10040 Leini, (Torino) - Italy
Tel. +39 011 997 9111
Fax. +39 011 997 9350
Toll-Free: 00 800 234 234 00
vpt-customer-care@agilent.com

Customer Support & Service
NORTH AMERICA:
Toll Free: 800 882 7426
vpl-ra@agilent.com
Lexington-service@agilent.com

EUROPE:
Toll-Free: 00 800 234 234 00
vpt-customer-care@agilent.com

PACIFIC RIM:
please visit our website for individual office
information
<http://www.agilent.com/chem/vacuum>

Worldwide Web Site, Catalog and Order
Online:
www.agilent.com/chem/vacuum

Representatives in most countries