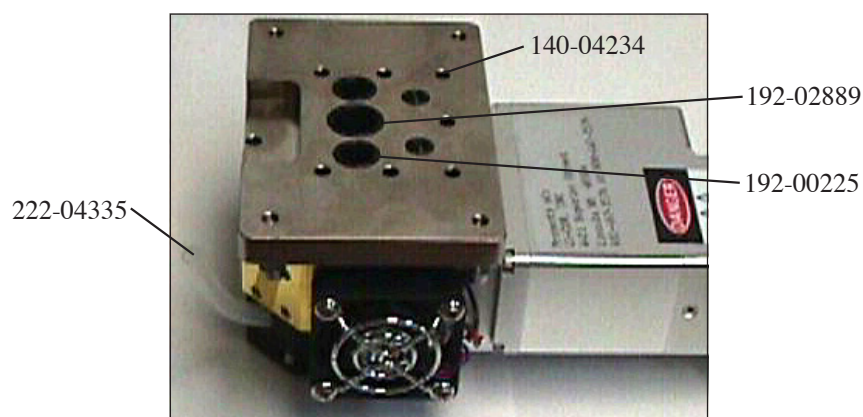
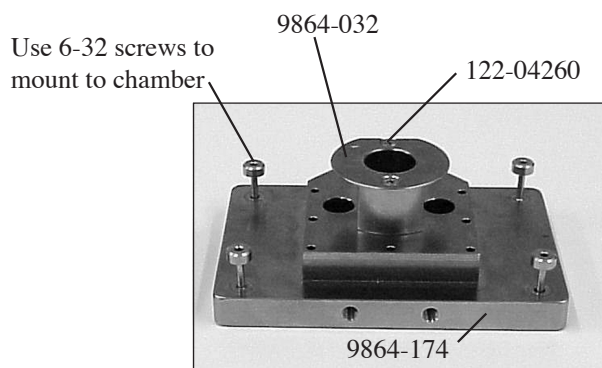


# 6400-19 Custom Chamber Kit

PPS-234

Quantity	Part Number	Description
2	192-02889	Large O-ring
4	192-00225	Small O-ring
4	122-04260	Chamber fan inlet screws
12	140-04234	Mounting plate hex screws
12 inches	222-04335	Exhaust/match tubing
1	309-04263	Thermocouple connector, male
1	434-09031	Thermocouple type E, .003" bead
3.0 feet	212-02314	Teflon tape
1	9964-080	Propafilm, 23 cm wide
1	9864-174	Mounting plate - dimensional drawing available <a href="#">here</a>
1	9864-032	Chamber fan inlet



---

### **LI-COR Biosciences - Global Headquarters**

4647 Superior Street  
Lincoln, Nebraska 68504  
Phone: +1-402-467-3576  
Toll free: 800-447-3576  
Fax: +1-402-467-2819  
envsales@licor.com • envsupport@licor.com • www.licor.com/env

### **Regional Offices**

#### **LI-COR GmbH, Germany**

*Serving Andorra, Albania, Cyprus, Estonia, France, Germany, Iceland, Latvia, Lithuania, Liechtenstein, Malta, Moldova, Monaco, San Marino, Ukraine, and Vatican City.*  
LI-COR Biosciences GmbH  
Siemensstraße 25A  
61352 Bad Homburg, Germany  
Phone: +49 (0) 6172 17 17 771  
Fax: +49 (0) 6172 17 17 799  
envsales-gmbh@licor.com • envsupport-gmbh@licor.com

#### **LI-COR Ltd., United Kingdom**

*Serving Denmark, Finland, Ireland, Norway, Sweden, and UK.*  
LI-COR Biosciences UK Ltd.  
St. John's Innovation Centre  
Cowley Road  
Cambridge  
CB4 0WS  
United Kingdom  
Phone: +44 (0) 1223 422102  
Fax: +44 (0) 1223 422105  
envsales-UK@licor.com • envsupport-UK@licor.com

LI-COR Distributor Network: [www.licor.com/env/distributors](http://www.licor.com/env/distributors)

