



Red rubber Z12,657-8  
White rubber Z16,736-3

This set is intended to provide the chemist with a full range of airtight closures for most laboratory glassware including  $\text{F}$  joints, storage bottles, test tubes, and round-bottom and Erlenmeyer flasks. Each septum is identified by the number appearing on the upper surface when the sleeve is pulled down over the lip of the stoppered container.

**Suba•Seal No./Size** ( $\text{F}$  size)

*Catalog Number*

Each size, along with the Aldrich catalog number, Suba•Seal number and container neck internal diameter, is represented in its location within the carrying case to facilitate reordering.

<b>17/9.5mm</b> <i>Z12,457-5</i>	<b>57/25.5mm</b> ( $\text{F}$ 29/42) <i>Z12,457-5</i>		<b>9/6.5mm</b> <i>Z12,455-9</i>
<b>53/24.0mm</b> <i>Z12,467-2</i>		<b>13/8.0mm</b> ( $\text{F}$ 10/30) <i>Z12,456-7</i>	<b>37/17.5mm</b> ( $\text{F}$ 19/22) <i>Z12,462-1</i>
<b>49/22.0mm</b> <i>Z12,466-4</i>		<b>45/20.5mm</b> ( $\text{F}$ 24/40) <i>Z12,465-6</i>	
<b>41/19.0mm</b> <i>Z12,464-8</i>		<b>33/16.0mm</b> <i>Z12,461-3</i>	
<b>21/21.0mm</b> <i>Z12,458-3</i>	<b>29/14.0mm</b> <i>Z12,460-5</i>		<b>25/12.5mm</b> ( $\text{F}$ 14/20) <i>Z12,459-1</i>

Questions ? or for ordering information call 800-558-9160 (USA/Canada) or 414-273-3850



**ALDRICH®**

*Organics and Inorganics for  
Chemical Synthesis*

1001 West Saint Paul Ave., Milwaukee, WI 53233

Telephone 414-273-3850 • 800-231-8327

Fax 414-273-4979 • 800-962-9591

Internet <http://www.sial.com/aldrich>

E-mail [aldrich@sial.com](mailto:aldrich@sial.com)

© 1997 Sigma-Aldrich Co. Printed in the USA. Aldrich is a member of the Sigma-Aldrich Family. Sigma-Aldrich warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product for its particular use. See reverse side of invoice or packing slip for additional terms and conditions of sale.