

Agilent Rack Adapter 3 Units

Product Name	Part Number
Rack Adapter 3 Units	9699191

The rack adapter shown in figure 1 is a 19" modular chassis system based on DIN 41494 part 5 manufactured by "Eurocard".

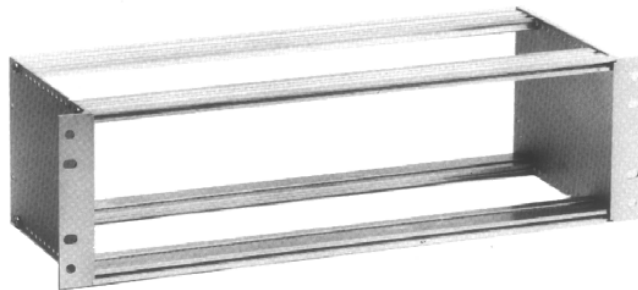


Figure 1 - Rack adapter 3 units

The modular chassis accepts standard 3U control units and provides mechanical mounting in cases or racks corresponding to the dimensional requirements of 19" front panels (DIN 41494 part 1).

The basic structure of this modular chassis system provides a further advantage.

The aluminum side panels are connected cross extrusions by four screws to give low-weight strength and stability.

This system reduces mounting times and cuts mechanical systems down to a minimum.

Low chassis height provides easy mounting on chassis runners.

Telescopic slides may be fixed to standard mounting holes on the side panels.

Instructions for use

Assemble the modular chassis as of figure 2 with the cross extrusion (3) so far enough to support the control units to be installed.

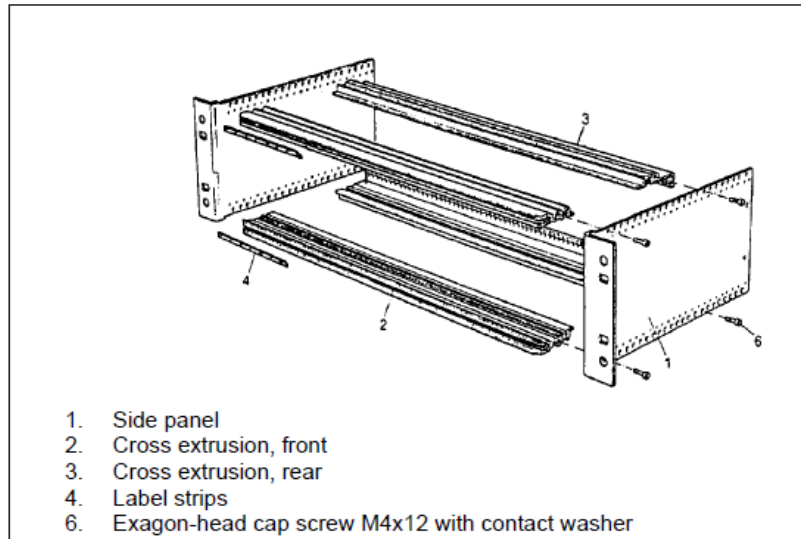


Figure 2 - Modular chassis assembly

The outline dimensions for the modular chassis are shown in figure 3.

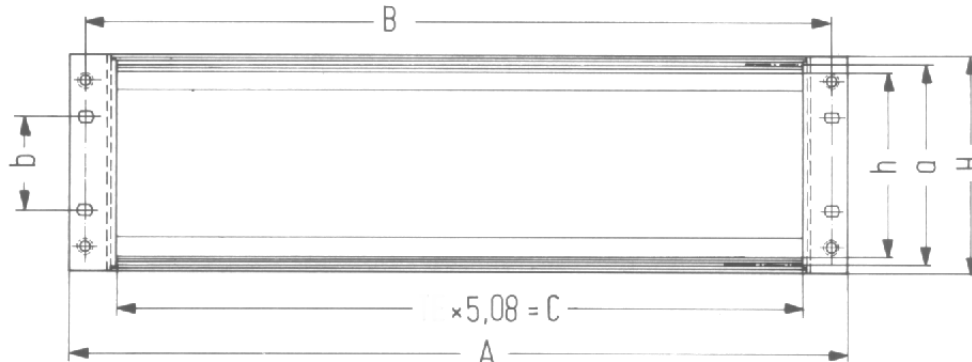


Figure 3 - Outline dimensions

Instructions for use

Mounting height 3U (dimensions mm/(inches))							
HP	A	B	C	H	a	b	h
84	<u>482.6</u> (19)	<u>465</u> (18.3)	<u>426.72</u> (16.8)	<u>132.5</u> (5.21)	<u>122.5</u> (4.82)	<u>57.15</u> (2.25)	<u>112.5</u> (4.42)

Where:

U (mounting height) is multiple of the height unit $1U = 44.45 \text{ mm} = 1\frac{3}{4}"$

HP (mounting width) is a multiple of the horizontal pitch (HP)

maximum of 84 HP

1 HP = 5.08 mm

Total width of the modular chassis = 482.6 mm = 19".

Instructions for use

Disposal

The following symbol is applied in accordance with the EC WEEE (Waste Electrical and Electronic Equipment) Directive. This symbol (valid only in countries of the European Community) indicates that the product it applies to must NOT be disposed of together with ordinary domestic or industrial waste but must be sent to a differentiated waste collection system. The end user is therefore invited to contact the supplier of the device, whether the Parent Company or a retailer, to initiate the collection and disposal process after checking the contractual terms and conditions of sale.



For more information refer to:

<http://www.agilent.com/environment/product/index.shtml>

Customer Support & Service:

NORTH AMERICA:

Toll Free: 800 882 7426

vp1-ra@agilent.com

Lexington-service@agilent.com

EUROPE:

Toll Free: 00 800 234 234 00

vp1-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

This information is subject to change without notice.

© Agilent Technologies, Inc. 2026

Printed in Italy, 03/2026

87-900-835-01 (C.00)