



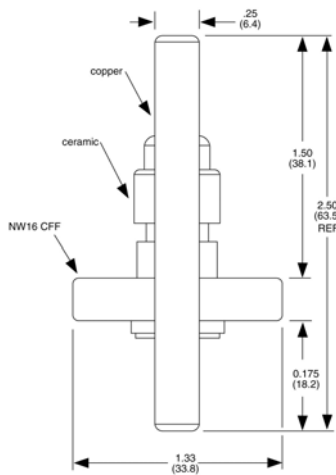
Agilent electrical feedthroughs provide reliable service in high and ultra high vacuum systems. Large, sturdy ceramic-to-metal seal construction provides maximum strenght and dependable performance over a wide range of vacuum applications.

All dimensions inches (mm)

75A feedthrough



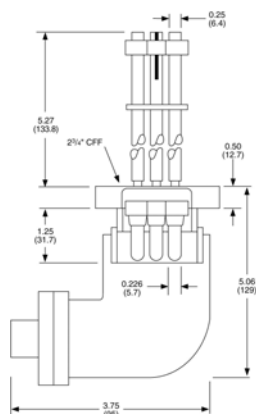
Model 954-5141



50A feedthrough



Model 954-5030



## SPECIFICATIONS

Feedthrough model number	954-5141	954-5030
Voltage rating (below 10 <sup>-4</sup> Torr/mbar)	2000V maximum	1kV pin to pin, max, 2kV pin to ground, max
Current rating	75A maximum	50A per pin maximum
Temperature bakeout range	-196°C to + 450°C	
Conductor	0.25" (6.4 mm) Ø OFHC copper rod	0.25" (6.4 mm) Ø nickel plated iron
Mounting flange	1.33" OD (NW 16) Mini ConFlat® flange	2.75" OD (NW35) ConFlat flange
Mating connectors vacuum external side	Push-on gold plated copper alloy, 150A, bakeable to +450°C, not insulated (Agilent 953-0074, not provided)	-----
Mating connector vacuum side	-----	Accepts wire or rod up to 0.102" OD (2.6 mm) (AWG 10)
Mating connector external side	-----	4 pins, 90° angle connector, provided (consists of: connector MS 3108B-32-17S Agilent 53-823217-00, clamp and bushing MS 3057-20A Agilent 51-969727-00)
Body material	Stainless steel	
Weight	1 lb - 0.45 kg	6 lb - 2.27. kg

## DISPOSAL

### Meaning of the "WEEE" logo found in labels

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.

