

**Instructions for use**

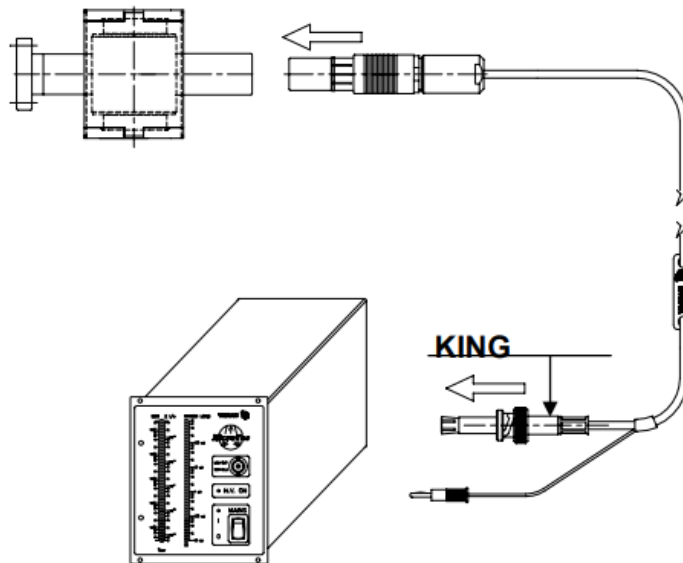
## Agilent H.V. Radiation Resistant Cable with Interlock/King

Product Name	Part Number
H.V. Radiation Resistant Cable with Interlock/King, 4 m	9290706



### SPECIFICATIONS

Operating voltage	7,5 kV.DC
Max temperature	-20°C / +220°C
Radiation resistance	10 E 9 rads
Minimum bending radius	60 mm
Net weight	lbs 0.95 (Kg 0.43)



## 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  
[vpl-ra@agilent.com](mailto:vpl-ra@agilent.com)  
[Lexington-service@agilent.com](mailto:Lexington-service@agilent.com)

##### EUROPE:

Toll Free: 00 800 234 234 00  
[vpt-customer-care@agilent.com](mailto: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](http://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

87-900-099-01 (C.00)