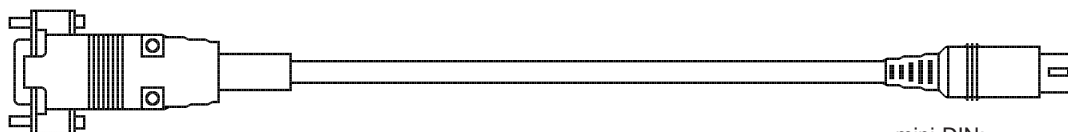


Model 2128 Fraction Collector Cables


Catalog Numbers
731-8286, 731-8290, 731-8291
731-8287, 731-8288

Econo System Cable #15, DB-15 to mini-DIN
Catalog Number 731-8286

Connections:

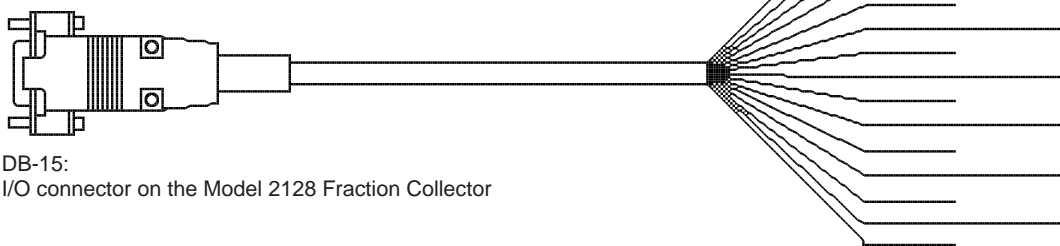


DB-15:
I/O connector on the Model 2128 Fraction Collector

mini-DIN:
 connector on the Econo System Controller

Model 2128 Accessory Cable, DB-15 to Bare Wires
Catalog Number 731-8290

Connections:



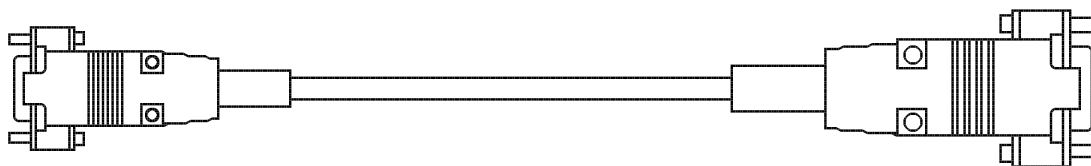
DB-15:
I/O connector on the Model 2128 Fraction Collector

1. Black
2. White
3. Red
4. Green
5. Orange
6. Blue
7. Wht/Blk
8. Red/Blk
9. Grn/Blk
10. Org/Blk
11. Blu/Blk
12. Blk/Wht
13. Red/Wht
14. Grn/Wht
15. Blu/Wht

Bare Wires:
Non Bio-Rad equipment

Model 2128 HPLC SI Cable, DB-15 to HPLC System Interface
Catalog Number 731-8291

Connections:

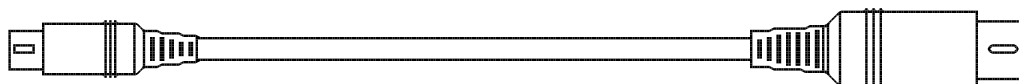


DB-15:
I/O connector on the Model 2128 Fraction Collector

HPLC:
Digital I/O connector on Bio-Rad's HPLC System Interface

Econo System Cable #16, Mini-DIN to DIN
Catalog Number 731-8287

Connections:



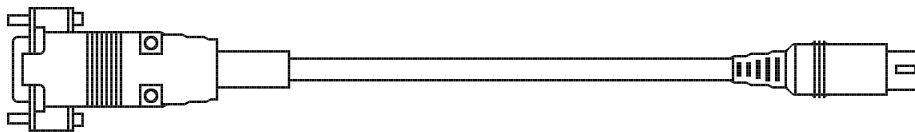
Mini-DIN:
REC connector on the Model 2128 Fraction Collector

DIN:
Channel 1 port on Bio-Rad Econo Recorder Models 1325,1326,1327

Model 2128 HPLC II Cable, 15 D to Bio-Rad's HPLC Instrument Interface

Catalog Number 731-8288

Connections:



15 D:
I/O connector on the Model 2128 Fraction Collector

Mini-DIN:
I/O connector on the Bio-Rad's HPLC Instrument Interface



**Bio-Rad
Laboratories**

*Life Science
Group*

Bio-Rad Laboratories Main Office, 2000 Alfred Nobel Drive, Hercules, California 94547, Ph. (510) 741-1000, Fx. (510) 741-1060 •
Eastern Regional Office, 85A Marcus Dr., Melville, New York 11747, Ph. (516) 756-2575, Fx. (516) 756-2594 •
Also in: North Ryde, Australia, Ph. 02-805-5000, Fx. 02-805-1920 • **Wien, Austria**, Ph. (1) 877 89 01, Fx. 0222-876 56 29 • **Nazareth, Belgium**,
Ph. 09-385 55 11, Fx. 09-385 65 54 • **Mississauga, Canada**, Ph. (905) 624-0713, Fx. (905) 624-3019 • **Beijing, China**, (01) Ph. 2563148, Fx. (01) 2564308 •
Ivry sur Seine Cedex, France, Ph. (1) 49 60 68 34, Fx. (1) 46 71 24 67 • **München, Germany**, Ph. 089 31884-0, Fx. 089 31884-100 • **Milano, Italy**,
Ph. 02-21609.1, Fx. 02-21609.399 • **Tokyo, Japan**, Ph. 03-3534-7515 Fx. 03-3534-8027 • **Veenendaal, The Netherlands**, Ph. 08385-40666,
Fx. 08385-42216 • **Auckland, New Zealand**, Ph. 09-443 3099, Fx. 09-443 3097 • **Kowloon, Hong Kong**, Ph. 7893300, Fx. 7891257 • **Flamingo Valley,
Singapore**, Ph. (65) 443 2529, Fx. (65) 442 1667 • **Solna, Sweden**, Phone 46 (0) 8 735 83 00, Fx 46 (0) 735 54 60 • **Madrid, Spain**, Ph. (91) 661 70 85,
Fx. (91) 661 96 98 • **Glattbrugg, Switzerland**, Ph. 01/810 16 77, Fx. 01/810 19 33 • **Hemel Hempstead, United Kingdom**, Ph. 0800 181134, Fx. 0442 259118

SIG 093093 Printed in USA

4006009 REV B