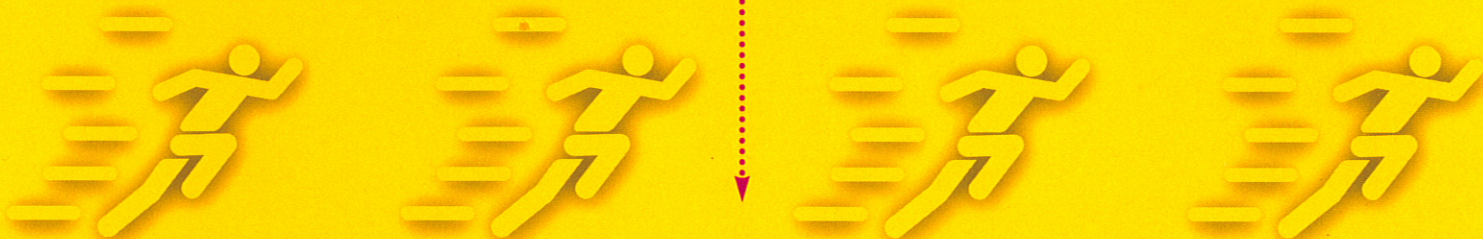


high performance

Chromatography Components

low pressure



Models:

EP-1 Econo Pump

EM-1 Econo UV Monitor

EG-1 Econo Gradient Monitor

2110 Fraction Collector

BIO-RAD