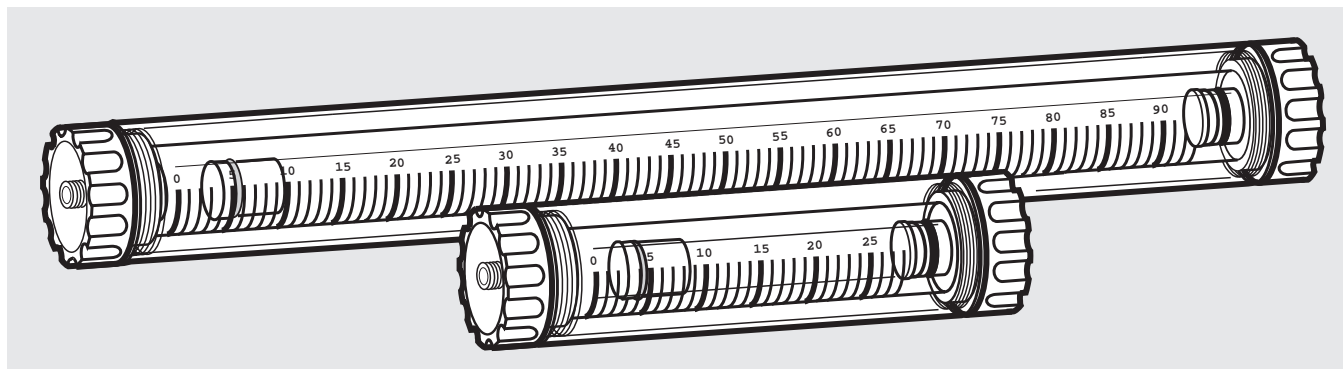


BioLogic DynaLoop 25 and 90 Sample Loops



The DynaLoop Sample Loop performs a wide range of sample injection volumes and options, making the BioLogic System more flexible than ever. The DynaLoop, available in 25 ml and 90 ml sizes, is used to load large volumes of sample directly onto chromatography columns. Injections can be made either manually or automatically, with the ease of simple fixed-loop injection.

The DynaLoop is a “dynamic” sample loop consisting of a high pressure glass tube (up to 1,000 psi), with an internal sliding seal that partitions sample from buffer. As the DynaLoop is loaded with sample, the sliding seal moves along as buffer is displaced. When sample is injected, the BioLogic gradient pump fills the DynaLoop with buffer, forcing the sliding seal in

the opposite direction, pushing the sample out of the loop and onto the column.

When plumbed directly to the AV7-3 or V7-3 BioLogic System injection valves, the DynaLoop functions just like a normal sample loop. Injection valve operation and sample loading are controlled automatically, simplifying the sample injection process and insuring precise sample loading and gradient formation.

The DynaLoop can be filled with sample manually with a syringe through valve port 5 or automatically using the model EP-1 Econo Pump (auxiliary pump, Figure 2).

All DynaLoop parts are user servicable which translates to inexpensive, easy-to-replace worn parts.

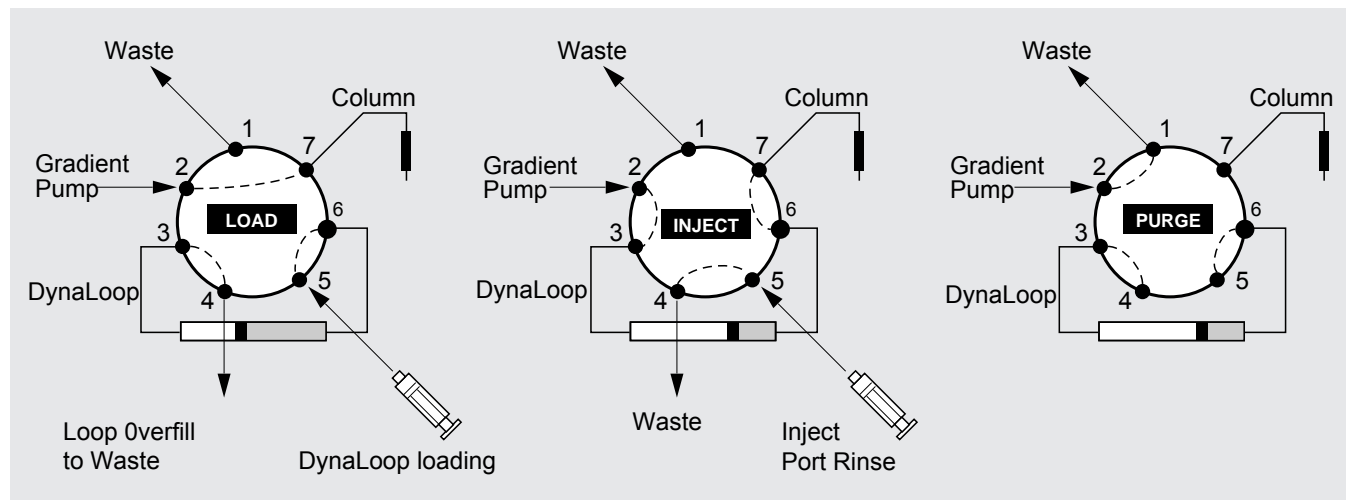


Fig. 1. AV7-3 or V7-3 Injection Valve plumbing with a DynaLoop.

DynaLoop Specification

Sizes	25 and 90 ml
Max. operating pressure	1,000 psi
Min. press. for seal movement	4 psi
Wetted materials	Tefzel, Ryton PTFE, PEEK, glass, Buna-N

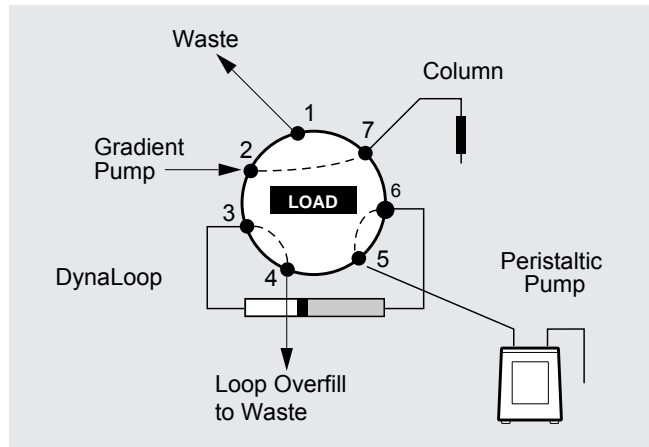


Fig. 2. Auxiliary pump tubing connections.

Ordering Information

Catalog Number	Product Description	Price
DynaLoops		
750-0451	DynaLoop Kit, includes a 25 ml DynaLoop, DynaLoop Parts Kit, and Instruction Manual	
750-0452	DynaLoop Kit, includes a 90 ml DynaLoop, DynaLoop Parts Kit, and Instruction Manual	
Spare Parts		
750-0450	DynaLoop Parts Kit, includes 4 end cap o-rings, 1 sliding seal o-ring, 1 filter, 4 collet/nut fittings, 4 1/8" ferrules, 5 1/4-28 nuts and ferrules, 1/8" PTFE tubing, 10 ft	

BIO-RAD

Bio-Rad
Laboratories

Life Science
Group

Bio-Rad Laboratories Main Office, 2000 Alfred Nobel Drive, Hercules, California 94547, Ph. (510) 741-1000, Fx. (510) 741-5800
 Also in: Reagents Park, Australia, Ph. 02-9914-2800, Fx. 02-9914-2888 Wien, Austria, Ph. (1) 877 89 01, Fx. (1) 876 56 29 Nazareth, Belgium, Ph. 09-385 55 11, Fx. 09-385 65 54
 Mississauga, Canada, Ph. (905) 712-2771, Fx. (905) 712-2990 Beijing, China, Ph. (01) 2046622, Fx. (01) 2051876 Copenhagen, Denmark, Ph. 39 17 9947, Fx. 39 27 1698
 Espoo, Finland, Ph. 90 804 2200, Fx. 90 804 1100 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
 New Delhi, India, Ph. 91-11-461-0103, Fx. 91-11-461-0765 Milano, Italy, Ph. 02-21609.1, Fx. 02-21609.399 Tokyo, Japan, Ph. 03-5811-6270 Fx. 03-5811-6272 Veenendaal,
 The Netherlands, Ph. 0318-540666, Fx. 0318-542216 Auckland, New Zealand, Ph. 09-443 3099, Fx. 09-443 3097 Kowloon, Hong Kong, Ph. 7893300, Fx. 7891257 Singapore,
 Ph. (65) 272-9877, Fx. (65) 273-4835 Solna, Sweden, Ph. 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. 01/809 55 55, Fx. 01/809 55 00 Hemel Hempstead, United Kingdom, Ph. 0800 181134, Fx. 01442 259118