

## Technical Bulletin

### AL-249

## Aldrich Sure/Pac Station Instructions

### Station for Non-Corrosive Liquefied Gases, Brass Regulator, Z566446

#### Specifications:

**Regulator:** Single stage, panel-mounted brass body with stainless steel pressure diaphragm, 1- $\frac{3}{4}$  inch diameter gauge, 300 psig maximum inlet pressure, 0 to 15 psig delivery pressure, PTFE-lined, braided stainless steel extension hose, brass swivel fitting and adapter for  $\frac{1}{4}$  inch NPTM Sure/Pac valve threads.

**Cabinet:** 10 in. W x 4 in. D x 3  $\frac{1}{2}$  in. H, Weight: 4 lb, without cylinders.



#### Components:

1. Station base
2. Brass regulator, includes brass outlet valve
3. PTFE-lined stainless hose
4. Grommet, rubber
5. Brass swivel fitting and adapter for  $\frac{1}{4}$  inch NPTM Sure/Pac valve threads

Sigma brand products are sold through Sigma-Aldrich, Inc.

Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply.

Please see reverse side of the invoice or packing slip.

## Start Up:

Place the 275 mL or 900 mL Sure/Pac cylinder into the cylinder slot in the station.\* Connect the ¼ inch NPT cylinder valve adapter to the Sure/Pac cylinder outlet valve using a wrench to tighten. Make sure to use PTFE tape on the threads for a proper leak proof seal. Thread the swivel nut from the connection hose onto the valve adapter and tighten with a wrench. (Never use PTFE tape on the JIC flare fitting. This fitting is designed to seal using the flare seat.) Make sure the regulator adjusting knob is turned out (counter clockwise) and the regulator outlet valve is closed (clockwise), so there will be no flow of gas when the cylinder is opened. Connect your use line to the regulator outlet valve, ⅜ inch NPTM, using PTFE tape as a sealant.

Tighten outlet connection with a wrench. Open the Sure/Pac cylinder valve slowly so that it is fully open. Check the cylinder valve and outlet connections for leaks using Snoop® or other compatible leak detector. Re-tighten or re-do if necessary. Slowly turn the regulator adjusting knob clockwise to increase delivery pressure to the desired pressure. Open regulator outlet valve to allow gas to flow.

\*Use the optional Cool/Heat Coil (see Accessories section) to warm the gas cylinder with circulating hot water to maintain gas flow during dispensing.

## Shut Down:

Turn off cylinder valve. Turn regulator adjusting knob clockwise until a small amount of delivery pressure shows on the delivery gauge. This allows all line pressure to be released from the regulator. Open outlet valve and allow residual pressure from the regulator and hose to drain from the system. Allow all gas pressure to drain from the gauge. Close outlet valve. Always drain gases safely into a proper hood or vent.



## Express Service™ Program (for US customers only):

**Regulator Repair** is offered as a service to all Sigma-Aldrich customers. We offer this convenience in order to expedite the repair of your equipment. We are able to repair regulation equipment subject to normal wear and tear such as replacement pressure gauges, seals, o-rings, diaphragm replacement, fittings, needle valves, springs, etc.

**Regulator Replacement** is another service offered to Sigma-Aldrich customers. If you have a regulator that has been sent to us but can not be repaired by our service center due to costs associated with the repair, damage due to corrosion or some other valid reason, we will offer you a new replacement regulator for your application at an affordable cost.

The procedure for this program is explained on our website at [sigma-aldrich.com/gases](http://sigma-aldrich.com/gases) where you will find complete instructions. For additional information, call our Technical Services Department at **800-231-8327**.

## Accessories:

### Cool/Heat Coil, Z566454

Copper coil warms gas cylinder to maintain gas flow during dispensing. Connects to hot tap water or recirculator using ¼ inch i.d. tubing. Coil fits both 275 mL and 900 mL, 3-inch diameter Sure/Pac cylinders.

