

Technical Bulletin

AL-151 Gas Equipment Configurations

The recommended regulators, control valves, lecture-bottle stations, and Sure/Pac stations for use with Aldrich gas products are specified on our web site, sigma-aldrich.com. Click the **Related Products** tab for a gas listing to view the recommended information.

Do not use Aldrich gas regulators on cylinders that exceed 3000 psig.

Configurations for Mildly Corrosive or High-Purity Gases



Regulator designated for mildly corrosive or high-purity gases with a T-purge assembly and bonnet vent tube.

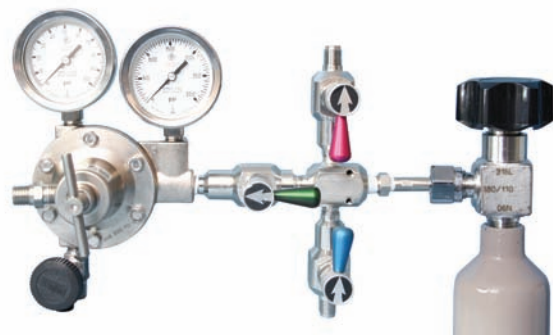


Control valve designated for mildly corrosive or high-purity gases to meter the flow of gas. **WARNING:** Gas-control valves do not control pressure in a closed system. Regulator or relief valves should be used to prevent the buildup of unsafe pressures that could lead to explosion.



Lecture-bottle station designated for mildly corrosive or high-purity gases.

Configurations for Extremely Corrosive Gases



Regulator with cross-purge assembly designated for extremely corrosive gases.

Configurations for Non-corrosive Gases



Regulator designated for non-corrosive gases with a check valve on the outlet side of the regulator to prevent back streaming of process gas.



Lecture-bottle station designated for extremely corrosive gases.



Lecture-bottle station designated for non-corrosive gases.



Control valve designated for extremely corrosive gases to meter the flow of gas. **WARNING:** Gas-control valves do not control pressure in a closed system. Regulator or relief valves should be used to prevent the buildup of unsafe pressures that could lead to explosion.



Control valve designated for non-corrosive gases to meter the flow of gas. **WARNING:** Gas-control valves do not control pressure in a closed system. Regulator or relief valves should be used to prevent the buildup of unsafe pressures that could lead to explosion.

Configurations for Liquefied Gases



Use the appropriate mini-regulator, brass or stainless steel



Use the appropriate hose barb adapter.



Sure/Pac station for non-corrosive liquefied gases