

AL-222

Technical Bulletin

Installation Notes for Pump N' Pour Adapter Cat. No. Z55,775-7

Because of variations in the design of the polyethylene bung gaskets that are used by the manufacturers of 20L solvent cans, the Pump N' Pour adapter may leak while pouring solvents.

To solve this problem, it is necessary to install the bung gasket that is supplied with the solvent can on the Pump N' Pour adapter.

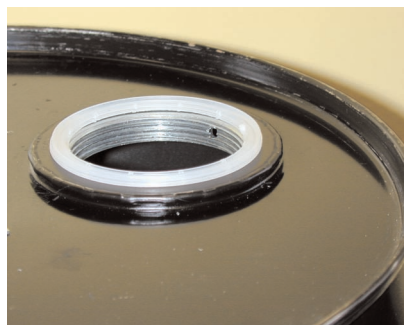
Follow these steps:

Remove the metal 2in. NPT bung from the 20L solvent can.

A bung wrench such as Aldrich **Cat. No. Z21,929-0** may be needed to loosen the bung initially.



Remove the flat polyethylene gasket from the bung.



Position the gasket on the threaded opening of the can.

Important Note: This gasket must rest in the recess of the threaded opening to seal properly. Verify that the gasket is positioned properly by moving the gasket side-side in the opening. You can feel that it is seated. If installed up-side down, you will not be able to seat the gasket in the threaded opening.

With the manufacturer's bung gasket in place, carefully thread the Pump N' Pour adapter into the opening and turn until hand tight. DO NOT use a wrench to tighten the adapter as this may strip the threads.



Observe that there is no leakage during the pouring operation.

Please contact the Technical Services department with any questions or problems regarding this Installation Note.

Telephone: 800-231-8327.