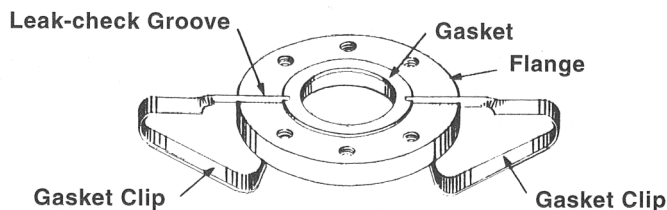


NOTICE: This document contains references to Varian. Please note that Varian, Inc. is now part of Agilent Technologies. For more information, go to [www.agilent.com/chem](http://www.agilent.com/chem).



**GASKET CLIP**  
MODEL NO. 953-5052  
P/N 5081-Q1455-001

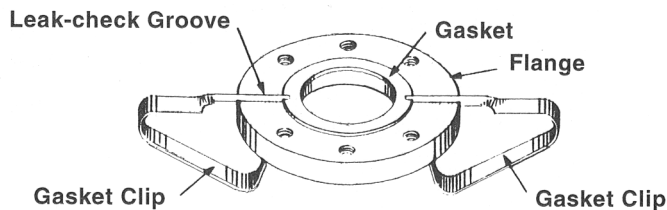


1. Use two gasket clips to hold a gasket in place during Conflat<sup>®</sup> flange assembly.
  2. Insert each clip in the flange's leak-check groove, as shown. The tip of the clip should not extend over the gasket more than 1/32 inch.
  3. Bolt flanges together finger-tight before removing the gasket clips.
- 04/2006 6999-05-030B

NOTICE: This document contains references to Varian. Please note that Varian, Inc. is now part of Agilent Technologies. For more information, go to [www.agilent.com/chem](http://www.agilent.com/chem).



**GASKET CLIP**  
MODEL NO. 953-5052  
P/N 5081-Q1455-001

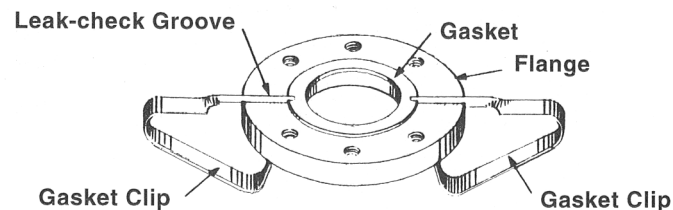


1. Use two gasket clips to hold a gasket in place during Conflat<sup>®</sup> flange assembly.
  2. Insert each clip in the flange's leak-check groove, as shown. The tip of the clip should not extend over the gasket more than 1/32 inch.
  3. Bolt flanges together finger-tight before removing the gasket clips.
- 04/2006 6999-05-030B

NOTICE: This document contains references to Varian. Please note that Varian, Inc. is now part of Agilent Technologies. For more information, go to [www.agilent.com/chem](http://www.agilent.com/chem).



**GASKET CLIP**  
MODEL NO. 953-5052  
P/N 5081-Q1455-001

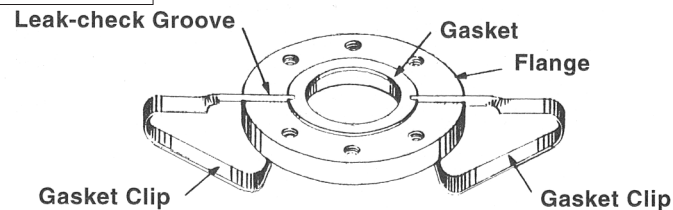


1. Use two gasket clips to hold a gasket in place during Conflat<sup>®</sup> flange assembly.
  2. Insert each clip in the flange's leak-check groove, as shown. The tip of the clip should not extend over the gasket more than 1/32 inch.
  3. Bolt flanges together finger-tight before removing the gasket clips.
- 04/2006 6999-05-030B

NOTICE: This document contains references to Varian. Please note that Varian, Inc. is now part of Agilent Technologies. For more information, go to [www.agilent.com/chem](http://www.agilent.com/chem).



**GASKET CLIP**  
MODEL NO. 953-5052  
P/N 5081-Q1455-001



1. Use two gasket clips to hold a gasket in place during Conflat<sup>®</sup> flange assembly.
  2. Insert each clip in the flange's leak-check groove, as shown. The tip of the clip should not extend over the gasket more than 1/32 inch.
  3. Bolt flanges together finger-tight before removing the gasket clips.
- 04/2006 6999-05-030B