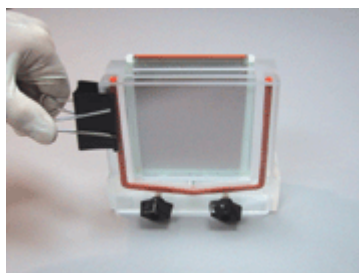


Thermo Scientific Owl P1-CST Multiple Gradient Gel Caster P1-CST Gel Casting Instructions

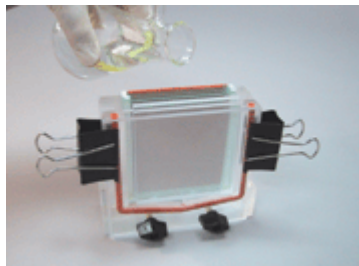
1. Assemble the gel cassette (s) and place into the caster.



2. Clamp the cassette in the caster.



3. Pour or gravity feed the acrylamide solution onto the cassette. Once the gel is polymerized, remove the cassette and place in a running device.



© 2012 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.

>> For more information, visit: www.thermoscientific.com/owlsci