













# ReadyPouch™ Agarose Gel

	<b>Package contents</b>	<b>Catalog Numbers</b> A25647 A25648	<b>Size</b> 10 × 100 mL 30 × 100 mL
	<b>Storage conditions</b>	Store at room temperature (do not freeze).	
	<b>Required materials</b>	<ul style="list-style-type: none"> <li>▪ 800–1200 W microwave oven</li> <li>▪ Agarose gel casting tray and combs</li> <li>▪ Molecular weight ladder</li> </ul>	
	<b>Timing</b>	<ul style="list-style-type: none"> <li>▪ Preparation: 90–130 seconds</li> <li>▪ Gel solidification: 20–40 minutes</li> </ul>	
	<b>Selection guide</b>	<a href="#">Agarose gel electrophoresis reagents</a> Go online to view related products.	
	<b>Product description</b>	<ul style="list-style-type: none"> <li>▪ Convenient agarose gel with TAE buffer in a pouch.</li> <li>▪ Ethidium bromide (0.625 mg/mL) is supplied in a plastic dropper. Use 2 drops/pouch for a final in-gel concentration of 0.5 µg/mL.</li> <li>▪ Can be used to separate DNA fragments ranging from 100 bp to 8 kb.</li> </ul>	
	<b>Important guidelines</b>	<ul style="list-style-type: none"> <li>▪ <b>CAUTION!</b> Contents will be hot after microwaving. Allow the pouch to cool briefly on the benchtop before pouring.</li> <li>▪ Reheating ReadyPouch™ Agarose Gels is not recommended. Buffer evaporation from heating multiple times can affect final agarose concentration.</li> <li>▪ Wear a lab coat, gloves, and eye protection.</li> <li>▪ Dispose of used ReadyPouch™ gel and ethidium bromide in accordance with the requirements of your EHS office and local regulations.</li> </ul>	
	<b>Online resources</b>	Visit our <a href="#">product page</a> for additional information and protocols. For support, visit <a href="http://www.lifetechnologies.com/support">www.lifetechnologies.com/support</a> .	



## Protocol

Steps		Procedure Details
1		<b>Vent</b> by tearing the pouch open along the dotted line.
2		<b>Heat</b> the pouch in a microwave oven on HIGH <ul style="list-style-type: none"> <li>▪ 90 seconds (&gt;1000 W oven) <b>or</b></li> <li>▪ 130 seconds (≤1000 W oven)</li> </ul>
3		<b>Add</b> 2 drops of ethidium bromide stain from the dropper supplied in the kit.
4		<b>Pour</b> the the agarose into a casting tray with the well comb in place.

### Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at [www.lifetechnologies.com/termsandconditions](http://www.lifetechnologies.com/termsandconditions). If you have any questions, please contact Life Technologies at [www.lifetechnologies.com/support](http://www.lifetechnologies.com/support).

### Important licensing information

This products may be covered by one or more Limited Use Label Licenses. By use of this product, you accept the terms and conditions of all applicable Limited Use Label Licenses.

### Disclaimer

LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

© 2014 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.