ABsolute qPCR Plate Seals

Catalog Number AB-1170

Pub. No. MAN0019011 Rev. B.0

Product description

The ABsolute qPCR Plate Seals is an optically clear, non-tacky pressure-sensitive adhesive developed specifically for qPCR and other fluorescence applications.

Contents and storage

Product	Cat. No.	Amount	Storage ^[1]
ABsolute qPCR Plate Seals	AB-1170	50 sheets	Keep cool and dryAvoid direct exposure to sunlight

^[1] It is recommended to use the seals within one year from date of purchase.

Product specifications

Sheet size	Length: 141.3 ± 1.6 mm (with tabs); 120.7 ± 1.6 mm (without tabs)		
	Width: 79.4 ± 1.6 mm		
Seal thickness	100 μm (without release liner)		
Temperature range	-80°C to 110°C		
Sealing	Bond strength is dependent on application pressure, so sufficient pressure must be applied. For optimal seal application, we recommend the use of the Adhesive Seal Applicator (Cat. No. AB-1391).		

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 at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.



Life Technologies Corporation | 5781 Van Allen Way | Carlsbad, CA 92008

For descriptions of symbols on product labels or product documents, go to thermofisher.com/symbols-definition.

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Important Licensing Information: This product 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.

©2022 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

