The Thermo Scientific ALPS 50 V microplate heat sealer is designed to provide sealing consistency in low-throughput environments. Heat sealing offers superior sample protection of storage and reaction plates in applications including compound storage, sample archiving and PCR.

Thermo Scientific ALPS™ 50 V Microplate Heat Sealer







Quality Sealing Options. With the semiautomated Thermo Scientific ALPS 50 V microplate heat sealer, secure tight seals around individual wells to eliminate sample loss through evaporation and cross contamination between wells. Choose from a wide range of heat seals, including piercable, optically clear and permanent, for any application.

Application Flexibility. Simple on-board touch pad controls adjust sealing temperature and time. Coupled with an audible alarm indicating sealing completion, the ALPS 50 V sealer provides optimal sealing control and ensures consistency.

Ergonomic Design. The counter lever handle design provides minimal strain when operating the instrument. Additionally, the ALPS 50 V sealer weighs only 7.2 kg (15.9 lbs) for ease of placement and relocation.

Knowledgeable Support. Our team has a strong reputation for supporting both equipment and consumables, and can demonstrate the ALPS 50 V sealer in your laboratory, as well as determine which seal is best for your application.



Thermo Scientific ALPS 50 V Microplate Heat Sealer

Dimensions (W x D x H)	220 cm x 321 cm x 425 cm	Applicable Standards	FCC Class A, CFR 47 Part 15 BS EN 61326:1998
Weight	15.9 lbs/ 7.2 kg	Power Supply	100-130 V AC, 50-60 Hz or 220-240 V AC, 50-60 Hz
Temperature Set Point Range	125 °C to 200 °C in 1 °C increments	Fuse	350 W
Time Set Point Range	1 to 9 seconds in 0.5 second increments	Power Rating	100-130 V - T3.15A (IEC 127) or 220-240V - T1.6A (IEC 127)
Operating Humidity Range	20-80% non-condensing	Mains Input Connector	IEC 320
Operating Temperature Range	18 °C to 30 °C		

Seals

Item No.	Description	Characteristics	Compatible Plate Material	Sealing Conditions	Seal Integrity Range	Applications	
AB-0559	Thermo-Seal	Foil laminate; very strong seal that can only be peeled using the Foil Stripper (see ordering information below); can be pierced with a single pipette tip or Pierce Plate (see ordering information below); good solvent resistance; recommended for use with DMSO	PP	1.5 to 2.5 seconds at 165 °C to 170 °C	-80 °C to 120 °C	Long term storage including storage at low temperatures; transportation and high temperature applications such as PCR, including use in waterbath thermal cyclers	
AB-0757	Easy Pierce	Foil material; easily pierced; peelable bond that can be removed by hand; second seal can be applied over existing seal; recommended for use with DMSO	PP, PE	0.5 to 3.0 seconds at 165 °C to 175 °C	-80 °C to 80 °C and up to 120 °C with cycler lid pressure	Well suited for 384-well formats where piercing is a useful method for sample retrieval; second seal can be applied over existing seal; suitable for PCR with screw-down lid	
AB-1720	Easy Pierce 20 µm	Thin foil material; easily pierced; peelable bond that can be removed by hand; good solvent resistance; recommended for use with DMSO	PP, PE	0.5 to 2.0 seconds at 165 °C to 175 °C	-80 °C to 80 °C and up to 120 °C with cycler lid pressure	Well suited for 384-well formats where piercing is a useful method for sample retrieval; suitable for PCR with screw-down lid	
AB-0745	Easy Peel	Foil laminate; difficult to pierce; peelable bond that can be easily removed by hand, but forms weld with PE plates; can be removed and new seal applied several times; some solvent resistance and suitable for DMSO use below 4 °C	PP, PE, COC	1.5 to 2.5 seconds at 165 °C to 170 °C	-200 °C to 90 °C and up to 120 °C with cycler pressure	Ideal for long term storage at low temperatures; seal can be applied, removed and a new seal applied several times; suitable for PCR with screw-down heated lid thermal cyclers only	
AB-1797	Clear Seal	Thin polymer material; able to be pierced; peelable seal when layered; good optical clarity; some solvent resistance	PP, PS, PE	1.5 seconds at 170 °C	-80 °C to 80 °C and up to 120 °C with cycler pressure	Suitable for fluorescence and colorimetric applications; suitable for PCR applications with screw-down or clip-down heated lid thermal cyclers only	
AB-0812	Clear Seal Diamond	Thin polymer material; peelable seal; good optical clarity; some solvent resistance	PP, PS, PE, COC	1.5 to 3.0 seconds at 170 °C	-80 °C to 120 °C	Ideal for fluorescence and colorimetric applications; suitable for PCR applications with screw-down or clip-down heated lid thermal cyclers only	
AB-0685	Clear Seal Strong	Clear laminate; weld seal that cannot be peeled or pierced; good optical clarity and solvent resistance; recommended for use with DMSO; clear for sample inspection	PP	1.5 to 2.0 seconds at 170 °C	-80 °C to 120 °C	Clear for sample inspection; due to strong unpeelable seal properties, it is suitable for storage and disposal of hazardous material; also suitable for PCR applications	

Ordering Guide

Item No.	Description	Qty	
AB-1443	ALPS 50 V Microplate Heat Sealer	Each	
AB-0563/1000	96-Well PCR Plate Carrier	Each	
AB-0724	384-Well PCR Plate Carrier	Each	
AB-0827	Standard Microplate Carrier	Each	
AB-0579	96-Pin Pierce Plate	Each	
AB-0592	Foil Stripper	Each	

© 2008 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.

