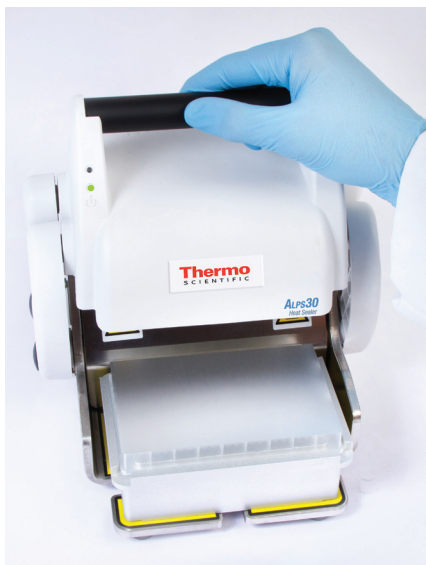


Thermo Scientific ALPS30 Manual Heat Sealer

Protect your precious samples simply and easily

The Thermo Scientific™ ALPS30 manual heat sealer is a hand-operated microplate heat sealer that allows you to protect your samples from evaporation and contamination when performing PCR or storing samples. Perfect for sealing small numbers of plates in low- to medium-throughput applications, the ALPS30 unit is small enough to keep on the bench and not take up a lot of space. Its simple, ergonomic manual operation provides quick, efficient sealing at your fingertips.



Sealing Flexibility.

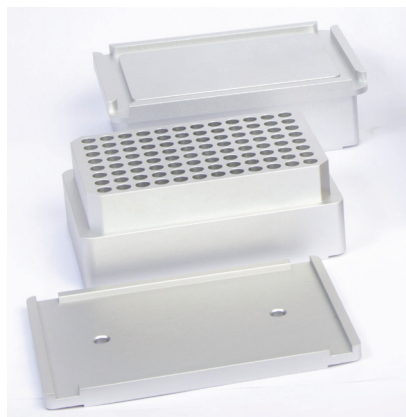
The ALPS30 sealer is compatible with virtually any height microplate or block, including PCR plates. The unit comes standard with three adapters to accommodate most plate types. The ALPS30 unit can be operated with the wide range of Thermo Scientific heat seals for any application, including pierceable, optically-clear and permanent seals.

Ultimate Protection.

The consistent heat seal applied by the ALPS30 sealer provides dependable protection to your PCR or sample storage plates. This seal prevents contamination and evaporation, affording you more reliable results.

Safe, Compact Design.

The ALPS30 sealer has a built-in thermostat to prevent overheating by triggering its automatic system shutoff. It features LED indicator lights showing when the power is on and when the unit is heating. The sealer unit is small and lightweight for easy carrying. Its small footprint means it doesn't take up a lot of space on your bench, so you can keep it handy for everyday use.



Thermo Scientific ALPS30 Manual Heat Sealer

Catalog No.	AB-0400
Dimensions (W x D x H), mm/in	220 x 250 x 230 / 8.7 x 9.8 x 9.1
Weight kg/lb	4 / 8.8
Operating Temperature Range	+18° to 30°C
Selecting Temperature Range	170°C (Non-adjustable) ±2°
Operating Humidity Range	20% to 80% non-condensing
Power Supply	110-120 / 220-230 VAC nominal IEC C13, 3-Core, 5A min
Fuse Ratings and Size	(115) T3.15AH 250V 20x5mm (230) T1.6AH 250V 20x5mm

Three adapters are included to accommodate most plate types:

- PCR Plates
- Deepwell plates and blocks
- Standard-height 96- or 384-format plates

Compatible Seals

Catalog 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 cyclers 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 cyclers 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 cyclers 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 cyclers 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

thermoscientific.com

© 2015 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.

Asia: Australia: 1300-735-292; New Zealand: 0800-933-966; China +86-21-6865-4588 or +86-10-8419-3588; China Toll-free: 800-810-5118 or 400-650-5118; Singapore +65-6872-9718; Japan: +81-3-5826-1616; Korea +82-2-2023-0640; Taiwan + 886-2-87516655; India: +91-22-6680-3000 **Europe:** Austria: +43-1-801-40-0; Belgium: +32-2-482-30-30; Denmark: +45-4631-2000; France: +33-2-2803-2180; Germany: +49-6184-90-6000; Germany Toll-free: 0800-1-536-376; Italy: +39-02-95059-554; Netherlands: +31-76-571-4440; Nordic/Baltic/CIS countries: +358-10-329-2200; Russia: +7-(812)-703-42-15; Spain/Portugal: +34-93-223-09-18; Switzerland: +41-44-454-12-12; UK/Ireland: +44-870-609-9203
North America: USA/Canada +1-585-586-8800; USA Toll-free: 800-625-4327
South America: USA sales support: +1-585-586-8800 **Countries not listed:** +49-6184-90-6000 or +33-2-2803-2000

Thermo
SCIENTIFIC

A Thermo Fisher Scientific Brand