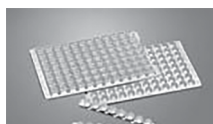




Sample storage and management

Storage plate sealer

Obtain a proper seal on plates, blocks and storage tube racks with the Thermo Scientific™ SuperSealer™ automatic sealer. Choose individual sealing options such as Thermo Scientific™ Matrix™ SepraSeal™ mats for liquid handling and single tube accessibility or seal entire plates and blocks for storage with Thermo Scientific™ Matrix™ CapMats™ storage plate cap mats. Effectively apply adhesive tape seals and ensure a tight seal with the included SuperSealer Adhesive Seal Mat.



Sealing simplicity

The SuperSealer storage plate sealer simplifies the sealing of plates, blocks and tube racks of virtually any height, with no adjustments needed, enabling a secure and consistent seal. Once the seal along with the plate, block or tube rack are placed into the drawer, simply close the door to initiate sealing.

Sealing flexibility

Designed to work with Matrix SepraSeal mats, CapMats and, with the SuperSealer adhesive seal mat, any pre-cut adhesive tape seal, you can be assured a proper seal regardless of sealing option.

Knowledgeable support

Our team has a strong reputation for supporting both equipment and consumables, and can demonstrate the SuperSealer plate sealer in your laboratory, detailing how it can address your labware sealing needs.

Product specifications	
Speed	10-15 seconds/plate, block or rack of tubes*
Dimensions (W x D x H)	6.5 "/16.5 cm x 12.5 "/31.75 cm x 13.5 "/34.3 cm
Weight	30 lbs/13.5 kg
Power supply	110-240 V

* The SuperSealer is designed to only work with full racks of tubes.



Ordering information

Description	Quantity	Cat. No.
Thermo Scientific SuperSealer		
SuperSealer	1 each	4110
SuperSealer adhesive seal mat	1 each	4111
Compatible Thermo Scientific matrix sealing options		
CapMat for 1 mL blocks, 96-well plates and snap racks	10/case	4411
CapMat for 1 mL blocks, 96-well plates and snap racks, sterile	10/case	4412
CapMat for 2 mL blocks, bulk	100/case	4420
CapMat for 2 mL blocks	10/case	4421
CapMat for 2 mL blocks, sterile	10/case	4422
CapMat for 500 µL, 0.75 ml and 1.4 mL storage tubes in latch racks, bulk	100/case	4430
CapMat for 500 µL, 0.75 ml and 1.4 mL storage tubes in latch rack	10/case	4431
CapMat for 500 µL, 0.75 mL and 1.4 mL storage tubes in latch racks, sterile	10/case	4432
SepraSeal mat	10 mats of 96 caps/case	4463
SepraSeal mats, sterile	10 mats of 96 caps/case	4464
SepraSeal mats, pre-split	10 mats of 96 caps/case	4465
SepraSeal mats, sterile, pre-split	10 mats of 96 caps/case	4466
SepraSeal cap removal tool	1 each	4469
Adhesive sealing tape for plates and block	100 seals/case	4417
Adhesive sealing tape for plates and blocks, dmso-resistant	100 seals/case	4418
Adhesive sealing tape for plates and blocks, foil	100 seals/case	4419

Learn more at thermofisher.com/supersealer

thermo scientific