



Thermo Scientific
Furnaces



Find a furnace
for your application

Thermo
SCIENTIFIC

Application	Recommended Furnaces	Heat Ranges	Furnace Type	Size	Embedded Heating Elements	Open Heating Elements
Alloying	Lindberg/Blue M heavy-duty 1200 °C box furnaces	🔥🔥🔥	Box	9-23 L		
	Lindberg/Blue M 1500 °C box furnace independent control	🔥🔥🔥🔥	Box	5.9 L	•	
	Lindberg/Blue M multi-purpose 1500 °C box furnaces	🔥🔥🔥🔥	Box	6 - 25 L	•	
	Lindberg/Blue M 1700 °C box furnaces, large chamber, integral control	🔥🔥🔥🔥🔥	Box	17 - 25.5 L		•
Analytical Processes	M104 muffle furnaces	🔥	Box	3.5 L	•	
	K114 chamber furnaces	🔥	Box	3.5 L	•	
Annealing	M104 muffle furnaces	🔥	Box	3.5 L	•	
	K114 chamber furnaces	🔥	Box	3.5 L	•	
	M110 muffle furnaces	🔥	Box	9 L	•	
	Lindberg/Blue M LGO 1200 °C box furnaces	🔥🔥🔥	Box	16.4 - 55.3 L	•	
	Lindberg/Blue M 1500 °C box furnace, independent control	🔥🔥🔥🔥	Box	5.9 L	•	
	Lindberg/Blue M multi-purpose 1500 °C box furnaces	🔥🔥🔥🔥	Box	6 - 25 L	•	
	Lindberg/Blue M 1200 °C split-hinge tube furnaces	🔥🔥🔥	Tube	30.5/61/91.4 cm (12/24/36 in.)		•
	Lindberg/Blue M 1500 °C general-purpose tube furnaces	🔥🔥🔥🔥	Tube	30.5 cm (12 in.)		•
	Lindberg/Blue M 1700 °C tube furnaces	🔥🔥🔥🔥🔥	Tube	30.5/61 cm (12/24 in.)		•
	Lindberg/Blue M 1200 °C crucible furnace, top loading	🔥🔥🔥	Crucible	5 in./7.5 in. ID, 8 in. Depth	•	
	Ashing	Thermolyne atmosphere controlled ashing furnaces	🔥	Box	14-45 L	•
Thermolyne premium large muffle furnaces		🔥🔥🔥	Box	14 L	•	
M104 muffle furnaces		🔥	Box	3.5 L	•	
K114 chamber furnaces		🔥	Box	3.5 L	•	
Lindberg/Blue M LGO 1200 °C box furnaces		🔥🔥🔥	Box	16.4 - 55.3 L	•	
Lindberg/Blue M heavy-duty 1200 °C box furnaces		🔥🔥🔥	Box	9-23L	•	
Lindberg/Blue M 1700 °C box furnaces, large chamber, integral control		🔥🔥🔥🔥🔥	Box	17 - 25.5 L		•
Lindberg/Blue M 1700 °C box furnaces independent control		🔥🔥🔥🔥🔥	Box	2.5-9 L		•
Atmosphere Processing	Thermolyne atmosphere controlled ashing furnaces	🔥	Box	14-45 L	•	
	Lindberg/Blue M 1500 °C general-purpose tube furnaces	🔥🔥🔥🔥	Tube	30.5 cm (12 in.)		•
	Lindberg/Blue M 1700 °C tube furnaces	🔥🔥🔥🔥🔥	Tube	30.5/61 cm (12/24 in.)		•
Baking	K114 chamber furnaces	🔥	Box	3.5 L	•	
	M110 muffle furnaces	🔥🔥	Box	9 L	•	
Bonding	Lindberg/Blue M 1700 °C box furnaces independent control	🔥🔥🔥🔥🔥	Box	2.5-9 L		•
Brazing	Lindberg/Blue M 1500 °C box furnace, independent control	🔥🔥🔥🔥	Box	5.9 L	•	
	Lindberg/Blue M Multipurpose 1500 °C box furnaces	🔥🔥🔥🔥	Box	6-25 L	•	
Calibration	Lindberg/Blue M mini-mite tube furnaces	🔥🔥	Tube	30.5cm (12 in.)		•
	Lindberg/Blue M 1200 °C split-hinge tube furnaces	🔥🔥🔥	Tube	30.5/61/91.4 cm (12/24/36 in.)		•
Ceramics	Lindberg/Blue M 1200 °C crucible furnace, top loading	🔥🔥🔥	Crucible	5 in./7.5 in. ID, 8 in. Depth	•	
Crystal Growing	Lindberg/Blue M 1200 °C split-hinge tube furnaces	🔥🔥🔥	Tube	30.5/61/91.4 cm (12/24/36 in.)		•
	Lindberg/Blue M 1700 °C tube furnaces	🔥🔥🔥🔥🔥	Tube	30.5/61 cm (12/24 in.)		•
Determination of Volatile and Suspended Solids	Thermolyne small benchtop muffle furnaces	🔥🔥	Box	1.3 - 2.1 L	•	
Drying	Lindberg/Blue M LGO 1200 °C box furnaces	🔥🔥🔥	Box	16.4 - 55.3 L	•	
Electronics	Lindberg/Blue M 1200 °C crucible furnace, top loading	🔥🔥🔥	Crucible	5 in./7.5 in. ID, 8 in. Depth	•	
Enameling	Lindberg/Blue M LGO 1200 °C box furnaces	🔥🔥🔥	Box	16.4 - 55.3 L	•	
Fusion	Lindberg/Blue M heavy-duty 1200 °C box furnaces	🔥🔥🔥	Box	9 - 23 L	•	
	Lindberg/Blue M 1700 °C box furnaces, large chamber, integral control	🔥🔥🔥🔥🔥	Box	17 - 25.5 L		•
Glass	Lindberg/Blue M 1200 °C crucible furnace, top loading	🔥🔥🔥	Crucible	5 in./7.5 in. ID, 8 in. Depth	•	
Gravimetric Analysis	Thermolyne small benchtop muffle furnaces	🔥🔥	Box	1.3 - 2.1 L	•	
	Thermolyne benchtop muffle furnaces	🔥🔥🔥	Box	2 - 5.8 L		•
	Thermolyne premium large muffle furnaces	🔥🔥🔥	Box	14 L	•	
	Thermolyne largest tabletop muffle furnaces	🔥	Box	45 L	•	
	Lindberg/Blue M 1100 °C tube furnaces	🔥🔥	Tube	61/91.4 cm (24/36 in.)		•

Application	Recommended Furnaces	Heat Ranges	Furnace Type	Size	Embedded Heating Elements	Open Heating Elements
Heat Treating	Thermolyne small benchtop muffle furnaces	🔥🔥	Box	1.3 - 2.1L	•	
	Thermolyne Industrial benchtop muffle furnaces	🔥🔥🔥	Box	2.2 L	•	
	Thermolyne benchtop muffle furnaces	🔥🔥🔥	Box	2 - 5.8 L		•
	Thermolyne largest tabletop muffle furnaces	🔥	Box	45 L	•	
	Lindberg/Blue M LGO 1200 °C box furnaces	🔥🔥🔥	Box	16.4 - 55.3 L	•	
	Lindberg/Blue M 1500 °C box furnace, independent control	🔥🔥🔥🔥	Box	5.9 L	•	
	Lindberg/Blue M Multipurpose 1500 °C box furnaces	🔥🔥🔥🔥	Box	6 - 25 L	•	
	Lindberg/Blue M 1100 °C tube furnaces	🔥🔥	Tube	61/91.4 cm (24/36 in.)		•
	Lindberg/Blue M 1200 °C split-hinge tube furnaces	🔥🔥🔥	Tube	30.5/61/91.4 cm (12/24/36 in.)		•
	Lindberg/Blue M 1500 °C general-purpose tube furnaces	🔥🔥🔥🔥	Tube	30.5 cm (12 in.)		•
Lindberg/Blue M 1700 °C tube furnaces	🔥🔥🔥🔥	Tube	30.5/61 cm (12/24 in.)		•	
Ignitions	Lindberg/Blue M heavy-duty 1200 °C box furnaces	🔥🔥🔥	Box	9 - 23L	•	
	Lindberg/Blue M 1700 °C box furnaces, Lindberg/Blue M large chamber, integral control	🔥🔥🔥🔥	Box	17 - 25.5 L		•
Incineration	K114 chamber furnaces	🔥	Box	3.5 L	•	
	M110 muffle furnaces	🔥🔥	Box	9 L	•	
Materials Testing	M104 muffle furnaces	🔥	Box	3.5 L	•	
Melting	Lindberg/Blue M LGO 1200 °C box furnaces	🔥🔥🔥	Box	16.4 - 55.3 L	•	
	Lindberg/Blue M 1500 °C box furnace, independent control	🔥🔥🔥🔥	Box	5.9 L	•	
	Lindberg/Blue M multi-purpose 1500 °C box furnaces	🔥🔥🔥🔥	Box	6 - 25 L	•	
	Lindberg/Blue M 1700 °C box furnaces independent control	🔥🔥🔥🔥	Box	2.5 - 9 L		•
	Lindberg/Blue M 1500 °C general-purpose tube furnaces	🔥🔥🔥🔥	Tube	30.5 cm (12 in.)		•
	Lindberg/Blue M 1200 °C crucible furnace, top loading	🔥🔥🔥	Crucible	5 in./7.5 in. ID, 8in. Depth	•	
Metallurgy	Lindberg/Blue M 1200 °C crucible furnace, top loading	🔥🔥🔥	Crucible	5 in./7.5 in. ID, 8in. Depth	•	
Metals and Ceramic Composites	Lindberg/Blue M 1700 °C box furnaces independent control	🔥🔥🔥🔥	Box	2.5 - 9L		•
Pyrolysis	Lindberg/Blue M mini-mite tube furnaces	🔥🔥	Tube	30.5 cm (12 in.)		•
Quantitative Analysis	Thermolyne benchtop muffle furnaces	🔥🔥🔥	Box	2 - 5.8L		•
	Thermolyne largest tabletop muffle furnaces	🔥	Box	45 L	•	
	Lindberg/Blue M 1100 °C tube furnaces	🔥🔥	Tube	61/91.4 cm (24/36 in.)		•
Sintering	Thermolyne benchtop muffle furnaces	🔥🔥🔥	Box	2 - 5.8 L		•
	Thermolyne largest tabletop muffle furnaces	🔥	Box	45 L	•	
	Lindberg/Blue M heavy-duty 1200 °C box furnaces	🔥🔥🔥	Box	9 - 23 L	•	
	Lindberg/Blue M 1700 °C box furnaces, large chamber, integral control	🔥🔥🔥🔥	Box	17 - 25.5 L		•
	Lindberg/Blue M 1700 °C box furnaces independent control	🔥🔥🔥🔥	Box	2.5 - 9 L		•
	Lindberg/Blue M mini-mite tube furnaces	🔥🔥	Tube	30.5 cm (12 in.)		•
	Lindberg/Blue M 1100 °C tube furnaces	🔥🔥	Tube	61/91.4 cm (24/36 in.)		•
	Lindberg/Blue M 1500 °C general-purpose tube furnaces	🔥🔥🔥🔥	Tube	30.5 cm (12 in.)		•
	Lindberg/Blue M 1700 °C tube furnaces	🔥🔥🔥🔥	Tube	30.5/61 cm (12/24 in.)		•
Superconductor Materials Research	Lindberg/Blue M 1200 °C crucible furnace, top loading	🔥🔥🔥	Crucible	5 in./7.5 in. ID, 8in. Depth	•	
Tempering	Lindberg/Blue M LGO 1200 °C box furnaces	🔥🔥🔥	Box	16.4 - 55.3 L	•	
Thermal Expansion	Lindberg/Blue M mini-mite tube furnaces	🔥🔥	Tube	30.5 cm (12 in.)		•
Viscosity Testing	Lindberg/Blue M mini-mite tube furnaces	🔥🔥	Tube	30.5 cm (12 in.)		•

🔥 100 °C - 1000 °C 🔥🔥 100 - 1100 °C 🔥🔥🔥 100 °C - 1200 °C 🔥🔥🔥🔥 500 - 1500 °C 🔥🔥🔥🔥🔥 500 - 1700 °C

Box furnaces Tube furnaces Crucible furnaces

thermofisher.com/furnaces

© 2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

Australia +61 39757 4300
Austria +43 1 801 40 0
Belgium +32 53 73 42 41
Chin.a +800 810 5118 or
+400 650 5118
France +33 2 2803 2180
Germany national toll free 0800 1 536 376
Germany in.ternational +49 6184 90 6000

In.dia toll free 1800 22 8374
In.dia +91 22 6716 2200
Italy +39 02 95059 552
Japan +81 3 5826 1616
Netherlands +31 76 579 55 55
New Zealand +64 9 980 6700
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 22
UK/Ireland +44 870 609 9203
USA/Canada +1 866 984 3766

Other Asian countries +852 2885 4613
Countries not listed +49 6184 90 6000

Thermo
SCIENTIFIC

A Thermo Fisher Scientific Brand