



Thermo Scientific
Furnaces



Find a furnace
by temperature range

Thermo
SCIENTIFIC

100 - 1000°C	🔥	K114 Chamber furnaces
		M104 muffle furnaces
		Thermolyne atmosphere controlled ashing furnaces
		Thermolyne largest tabletop muffle furnaces
100 - 1100°C	🔥🔥	M110 muffle furnaces
		Thermolyne small benchtop muffle furnaces
		Lindberg/Blue M mini-mite tube furnaces
		Lindberg/Blue M 1100°C tube furnaces
100 - 1200°C	🔥🔥🔥	Thermolyne premium large muffle furnaces
		Thermolyne benchtop muffle furnaces
		Thermolyne industrial benchtop muffle furnaces
		Lindberg/Blue M heavy-duty 1200°C box furnaces
		Lindberg/Blue M 1200°C split-hinge tube furnaces
500-1500°C	🔥🔥🔥🔥	Lindberg/Blue M 1200°C Crucible furnace, top loading
		Lindberg/Blue M multi-purpose 1500°C box furnaces
		Lindberg/Blue M 1500°C box furnaces, independent control
500-1700°C	🔥🔥🔥🔥	Lindberg/Blue M 1500°C general-purpose tube furnaces
		Lindberg/Blue M 1700°C box furnaces, large chamber, integral control
		Lindberg/Blue M 1700°C box furnaces, independent control
		Lindberg/Blue M 1700°C tube furnaces

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