

# Discover a furnace for your application

Furnaces in laboratory and industrial settings are used daily for a wide variety of applications. Thermo Scientific™ furnaces are designed to provide reliability, safety and ease of use. With a wide array of features, our versatile offering can help you find a solution to meet your needs. Choose from a variety of temperature control options, heating elements and furnace platforms to meet your analytical requirements. With Thermo Scientific furnaces, achieve consistent results you can depend on.



Application	Recommended furnaces	Heat range	Furnace type	Capacity range	Heating element	
					Embedded	Open
Ashing	Thermolyne Small Benchtop Muffle Furnaces	🔥🔥	Box	1 - 2 L	✓	
	Thermolyne Benchtop Muffle Furnaces	🔥🔥🔥	Box	2 - 5 L		✓
	Thermolyne Premium Large Muffle Furnaces	🔥🔥🔥	Box	14 L	✓	
	Thermolyne Largest Tabletop Muffle Furnaces	🔥🔥🔥	Box	45 L	✓	
	Lindberg/Blue M Moldatherm Box Furnaces	🔥🔥	Box	2 - 45 L	✓	
	Lindberg/Blue M LGO Box Furnaces	🔥🔥🔥	Box	16 - 55 L	✓	
Calibration and crystal growing	Lindberg/Blue M Mini-Mite Tube Furnaces	🔥🔥	Tube	12 in		✓
	Lindberg/Blue M 1200 °C Split-Hinge Tube Furnaces	🔥🔥🔥	Tube	12   24   36 in		✓
Determination of volatile and suspended solids	Thermolyne Small Benchtop Muffle Furnaces	🔥🔥	Box	1 - 2 L	✓	
	Thermolyne Benchtop Muffle Furnaces	🔥🔥🔥	Box	2 - 5 L		✓
	Lindberg/Blue M Moldatherm Box Furnaces	🔥🔥	Box	2 - 45 L	✓	
Drying, enameling, and tempering	Lindberg/Blue M Moldatherm Box Furnaces	🔥🔥	Box	2 - 45 L	✓	
	Lindberg/Blue M LGO Box Furnaces	🔥🔥🔥	Box	16 - 55 L	✓	
Gravimetric and quantitative analysis	Thermolyne Small Benchtop Muffle Furnaces	🔥🔥	Box	1 - 2 L	✓	
	Thermolyne Benchtop Muffle Furnaces	🔥🔥🔥	Box	2 - 5 L		✓
	Thermolyne Premium Large Muffle Furnaces	🔥🔥🔥	Box	14 L	✓	
	Thermolyne Largest Tabletop Muffle Furnaces	🔥🔥🔥	Box	45 L	✓	
	Lindberg/Blue M Moldatherm Box Furnaces	🔥🔥	Box	2 - 45 L	✓	
	Lindberg/Blue M LGO Box Furnaces	🔥🔥🔥	Box	16 - 55 L	✓	
	Lindberg/Blue M Mini-Mite Tube Furnaces	🔥🔥	Tube	12 in		✓
Heat treatment, annealing, sintering, and bonding	Lindberg/Blue M 1200 °C Split-Hinge Tube Furnaces	🔥🔥🔥	Tube	12   24   36 in		✓
	Thermolyne Small Benchtop Muffle Furnaces	🔥🔥	Box	1 - 2 L	✓	
	Thermolyne Benchtop Muffle Furnaces	🔥🔥🔥	Box	2 - 5 L		✓
	Thermolyne Premium Large Muffle Furnaces	🔥🔥🔥	Box	14 L	✓	
	Thermolyne Largest Tabletop Muffle Furnaces	🔥🔥🔥	Box	45 L	✓	
	Lindberg/Blue M Moldatherm Box Furnaces	🔥🔥	Box	2 - 45 L	✓	
	Lindberg/Blue M LGO Box Furnaces	🔥🔥🔥	Box	16 - 55 L	✓	
	Lindberg/Blue M Mini-Mite Tube Furnaces	🔥🔥	Tube	12 in		✓
Pyrolysis	Lindberg/Blue M 1200 °C Split-Hinge Tube Furnaces	🔥🔥🔥	Tube	12   24   36 in		✓
	Lindberg/Blue M Mini-Mite Tube Furnaces	🔥🔥	Tube	12 in		✓
Thermal expansion and viscosity testing	Lindberg/Blue M Mini-Mite Tube Furnaces	🔥🔥	Tube	12 in		✓
	Lindberg/Blue M 1200 °C Split-Hinge Tube Furnaces	🔥🔥🔥	Tube	12   24   36 in		✓

🔥 100 °C - 1000 °C   🔥🔥 100 - 1100 °C   🔥🔥🔥 100 °C - 1200 °C

Learn more at [thermofisher.com/furnaces](https://thermofisher.com/furnaces)

**thermo**scientific

**For Laboratory Use. It is the customer's responsibility to ensure that the performance of the product is suitable for customers' specific use or application.** © 2023 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. **EXT4649 0323**