



## Thermo Scientific X-ray Sources

### The power, resolution, and performance to fulfill a variety of applications

X-ray sources are deployed in a wide range of applications, for the testing, imaging, and inspection of items ranging from batteries, printed circuits boards and semi-conductor devices to animal fluoroscopy and dental prosthetics quality control. X-ray sources have the short wavelengths and high energies required to penetrate through many different materials and can be powered on and off at will, eliminating any potential for harm when not in use.

The portfolio of Thermo Scientific™ X-ray Sources includes 4 distinct model families, with multiple variants offering the flexibility to closely match source performance to specific requirements. Key determinants of choice include maximum voltage and minimum spot size. We offer sources with the ability to produce minimum spot sizes in the microfocus range: 1  $\mu\text{m}$  to 100  $\mu\text{m}$ .

## The right source for each application

Our portfolio of x-ray sources covers a wide range of specifications to help you identify the best source for your application.

Product family	Model	Tube	Max power	Max voltage	Max current	Min spot size	FOD	X-ray beam angle
PXS5	PXS5-822	822	8 W	80 kV	0.178 mA	15 µm	12.5 +/- 0.5 mm	34°
	PXS5-925	925	8 W	90 kV	0.180 mA	5 µm	12.0 ± 0.5 mm	40°
	PXS5-925-LV	925	8 W	70 kV	0.180 mA	5 µm	12.0 ± 0.5 mm	40°
	PXS5-928-AN	928	8 W	90kV	0.160 mA	4.5 µm	6.0 mm	108°
	PXS5-928	928	8 W	90 kV	0.160 mA	4.5 µm	6.0 mm	108°
	PXS5-928-LV03	928	3 W	60 kV	0.075 mA	<5 µm	6.0 mm	108°
PXS10	PXS10-16W	131	16 W	130 kV	0.320 mA	6 µm	14.0 ± 0.5 mm	53°
	PXS10-40W	131	40 W	130 kV	0.500 mA	6 µm	14.0 ± 0.5 mm	53°
	PXS10-65W	131	65 W	130 kV	0.500 mA	6 µm	14.0 ± 0.5 mm	53°
	PXS10-WB-16W	134	16 W	130 kV	0.320 mA	7 µm	10.0 ± 0.5 mm	115°
	PXS10-WB-40W	134	40 W	130 kV	0.500 mA	7 µm	10.0 ± 0.5 mm	115°
	PXS10-WB-65W	134	65 W	130 kV	0.500 mA	7 µm	10.0 ± 0.5 mm	115°
	PXS10-WBE-16W	134	16 W	130 kV	0.320 mA	7 µm	10.0 ± 0.5 mm	115°
	PXS10-WBE-40W	134	40 W	130 kV	0.500 mA	7 µm	10.0 ± 0.5 mm	115°
	PXS10-WBE-65W	134	65 W	130 kV	0.500 mA	7 µm	10.0 ± 0.5 mm	115°
PXS11	PXS11-100-35-Rohs	7010	7.5 W	75 kV	0.100 mA	45 µm	8.9 mm	34° nominal
PXS15	PXS15-NB	110kV-NB	25 W	110 kV	0.250 mA	5 µm	7.5 ± 1.0 mm	90°
	PXS15-WB	110kV-WB	25 W	110 kV	0.250 mA	7 µm	7.5 ± 1.0 mm	125°

\* All PXS10 models available in standard or lightweight (LW) shielding options



Thermo Scientific™ PXS5  
Microfocus X-Ray Source



Thermo Scientific™ PXS10  
Microfocus X-Ray Source



Thermo Scientific™ PXS11  
Microfocus X-Ray Source



Thermo Scientific™ PXS15  
Microfocus X-Ray Source

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

**thermo** scientific