



# PANNORAMIC MIDI | Versatile Whole Slide Imaging System

## *Multi-modal whole slide scanner for fixed or wet specimens!*

The fully-automated Pannoramic MIDI™ generates whole slide images in transmitted light or fluorescence, both modalities available in the same unit. Gather data in up to 9 different fluorescence channels quickly and at superior resolution. The functional slide loading design accommodates both fixed or wet specimens. The 12-slide capacity makes it a perfect fit for small to medium-sized laboratories.

All 3DHitech scanners feature industry-leading image resolution, speed and remote access capabilities. The Pannoramic Viewer™ software that comes with the system allows for efficient image viewing, annotation and archiving capabilities. More advanced image analysis can be performed with plug-in tools, including tools for cytoplasmic, membrane or nuclear measurements. If automated image analysis is required, Caliper's inForm™ advanced image analysis software provides an easy-to-use interface that relies on a robust learn-by-example method.

Share and access images via your own intranet, using the powerful and flexible CaseCenter™. Unlike other similar packages, it allows for easy customization, access to remotely gathered data, open teleconsultation for second opinions, and it can run invisibly behind your organization's LIS portal.

## Benefits

- Whole slide imaging
- Teleconsultation
- Education
- Archiving
- Image Analysis

## MIDI Features

### SLIDE LOADING

Ergonomic and flexible, the Panoramic MIDI automatically scans up to 12 slides in one run. The unique horizontal holder allows for both fixed and wet specimens. The system automatically detects the sample and digitizes it with high quality.

### FLUORESCENCE

The Panoramic MIDI supports fluorescence scanning (optional) accommodating up to 9 different fluorescence channels.

### SPEED AND VERSATILITY

The Panoramic MIDI can scan a 10x10mm area in less than a minute. It is equipped with an external scan camera that can easily be changed to adjust for the application, allowing for improvement in speed and/or quality.

### QUALITY

The Panoramic MIDI supports the best Carl Zeiss™ objectives, ranging from 20x to 40x magnification. We achieve superior pixel resolution of 0.23  $\mu\text{m}/\text{pixel}$  with a 20x objective.

### SOFTWARE

Caliper and 3DHistech's software product portfolio is one of the most comprehensive. Whether you are scanning tissue microarrays or conventional IHC, or you require semi-automated or fully automated quantitative measurements, or teaching applications, there is a solution to fit your application.

## Pannoramic Scanner Specifications

	DESK	MIDI	SCAN	P250
Slide Capacity	1	12	150	250
Slide Loading	Manual	Automatic	Automatic	Automatic
Available Resolution Range (and Equivalent Objective Magnification)	0.50 $\mu\text{m}/\text{pixel}$ (equiv 20x mag) 0.23 $\mu\text{m}/\text{pixel}$ (equiv 45x mag)	0.50 $\mu\text{m}/\text{pixel}$ (equiv 20x mag) 0.23 $\mu\text{m}/\text{pixel}$ (equiv 45x mag)	0.50 $\mu\text{m}/\text{pixel}$ (equiv 20x mag) 0.23 $\mu\text{m}/\text{pixel}$ (equiv 45x mag)	0.35 $\mu\text{m}/\text{pixel}$ (equiv 29x mag) 0.22 $\mu\text{m}/\text{pixel}$ (equiv 45x mag)
Scanning Time (15 X 15 Mm)	1 min 49 sec	1 min 49 sec	1 min 49 sec	54 sec
Fluorescence Modality	No	Yes	Yes	Yes
Slide Type	FIXED or WET	FIXED or WET	FIXED	FIXED
Dimensions (W X H X D)	10.6 x 9.8 x 11.8" (27 x 25 x 30 cm)	18.9 x 27.6 x 19.7" (48 x 70 x 50 cm)	24 x 35.4 x 27.2" (61 x 90 x 69 cm)	24 x 35.4 x 27.2" (61 x 90 x 69 cm)
Weight	24 lbs	51 lbs	57 lbs	101.2 lbs
Objective(S)	20x/0.8 NA 40x/0.95 NA	20x/0.8 NA 40x/0.95 NA	20x/0.8 NA 40x/0.95 NA	20x/0.8 NA 40x/0.95 NA