

Conclusion

The limiting resolution of the Quantum GX and Quantum FX imaging systems is market leading. The theoretical limiting resolution of 9 μm in the Quantum GX and ability to visually identify a 15 μm line in a bar pattern phantom is better than any other system currently available. This resolution is quite sufficient relative to the common needs for small animal imaging applications. Small anatomical structures such as trabecular bone and small vessels can be easily resolved with the resolutions achieved on either imaging system.^{12, 13}

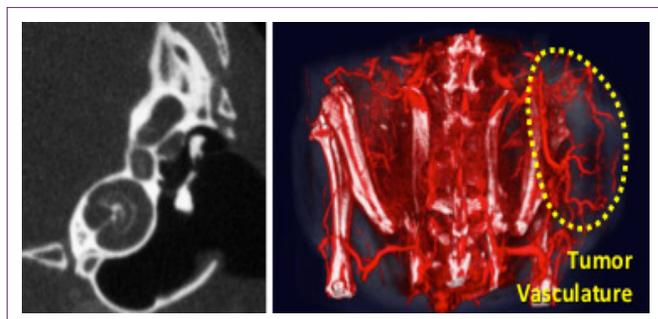


Figure 6: Images of (A) mouse inner ear structure and (B) tumor vasculature in a murine tumor model obtained using the Quantum GX.

References

1. J. F. Barrett and N. Keat, "Artifacts in CT: Recognition and Avoidance," *RadioGraphics*, pp. 1679–1691, 24AD.
2. J. a. Seibert and J. M. Boone, "X-ray imaging physics for nuclear medicine technologists. Part 2: X-ray interactions and image formation," *J. Nucl. Med. Technol.*, vol. 33, no. 1, pp. 3–18, 2005.
3. L. Y. Du, J. Umoh, H. N. Nikolov, S. I. Pollmann, T.-Y. Lee, and D. W. Holdsworth, "A quality assurance phantom for the performance evaluation of volumetric micro-CT systems," *Phys. Med. Biol.*, vol. 52, no. 23, pp. 7087–7108, 2007.
4. L. Arentsen and S. Hui, "Characterization of rotating gantry micro-CT configuration for the in vivo evaluation of murine trabecular bone," *Microsc. Microanal.*, vol. 19, pp. 907–13, 2013.
5. W. Stiller, M. Kobayashi, K. Koike, U. Stampfl, G. M. Richter, W. Semmler, and F. Kiessling, "Initial Experience with a Novel Low-Dose Micro-CT System," *Rofo*, vol. 179, no. 7, pp. 669–675, 2007.
6. S. N. Friedman, G. S. K. Fung, J. H. Siewerdsen, and B. M. W. Tsui, "A simple approach to measure computed tomography (CT) modulation transfer function (MTF) and noise-power spectrum (NPS) using the American College of Radiology (ACR) accreditation phantom," *Med. Phys.*, vol. 40, no. 5, p. 051907, 2013.
7. J. Rueckel, M. Stockmar, F. Pfeiffer, and J. Herzen, "Spatial resolution characterization of a X-ray microCT system," *Appl. Radiat. Isot.*, vol. 94, pp. 230–234, 2014.
8. O. Langner, M. Karolczak, G. Rattmann, and W. A. Kalender, "Bar and Point Test Patterns Generated by Dry-Etching for Measurement of High Spatial Resolution Micro-CT," in *IFMBE Proceedings, World Congress on Medical Physics and Biomedical Engineering*, 2009.
9. S. W. Smith, "The Scientist and Engineer's Guide to Digital Signal Processing," in *The Scientist and Engineer's Guide to Digital Signal Processing*, 1997, pp. 423–450.
10. J. P. Moy, "Signal-to-noise ratio and spatial resolution in x-ray electronic imagers: is the MTF a relevant parameter?," *Med. Phys.*, vol. 27, no. 1, pp. 86–93, 2000.
11. M. U. Ghani, Z. Zhou, L. Ren, M. Wong, Y. Li, B. Zheng, K. Yang, and H. Liu, "Investigation of spatial resolution characteristics of an in vivo microcomputed tomography system," *Nucl. Instruments Methods Phys. Res. Sect. A Accel. Spectrometers, Detect. Assoc. Equip.*, vol. 807, pp. 129–136, 2016.
12. J.-C. Tseng, A. Behrooz, J. Meganck, J. Kempner, and J. Peterson, "Using Quantum GX mCT system to investigate bone erosion and deformation in osteoarthritis rats" *PerkinElmer applications notes*, 2016.
13. N. M. Das, S. Hatsell, K. Nannuru, L. Huang, X. Wen, L. Wang, L.-H. Wang, V. Idone, J. A. Meganck, A. Murphy, A. Economides, and L. Xie, "In Vivo Quantitative Microcomputed Tomographic Analysis of Vasculature and Organs in a Normal and Diseased Mouse Model," *PLoS One*, vol. 11, no. 2, p. e0150085, Feb. 2016.