

Which Imaging Modality Is Best For Your Research?

ONCOLOGY

LIVER DISEASE

KIDNEY DISEASE

CARDIAC/CVD DISEASE

REPRODUCTIVE HEALTH

MSK/ADIPOSE

LUNG DISEASE



3D Tumor Volumes
Angiogenesis



Fibrosis.
Cancer, Fat



Vascularity and
Obstruction

Volume and
Stiffness/
Fibrosis



Cardiac Function

Vascular
Imaging



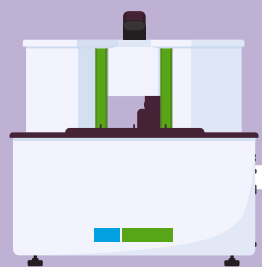
Pregnancy



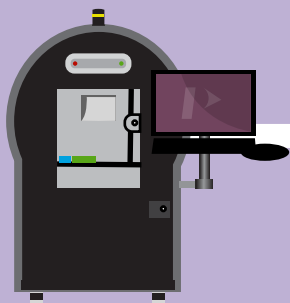
Muscle
Volume/
Fibrosis



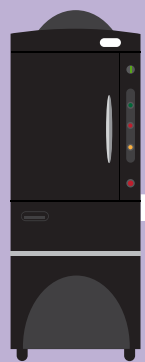
Volume,
Function
(Fibrosis
Cancer etc.)



Vega® Ultrasound

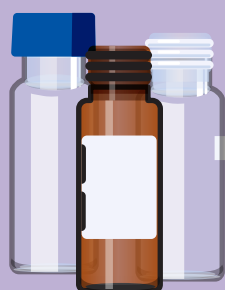


Quantum GX2 MicroCT



IVIS® Platform
Bioluminescence/Fluorescence

Inflammation, Gene activation, Cell Tracking, Biodistribution, Metastases, Metabolism, Disease Burden



Imaging Reagents
IVISbrite™, IVISense™, VesselVue®

For more information visit our website: www.perkinelmer.com/invivo or contact your sales representative.

Copyright ©2022, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

892808