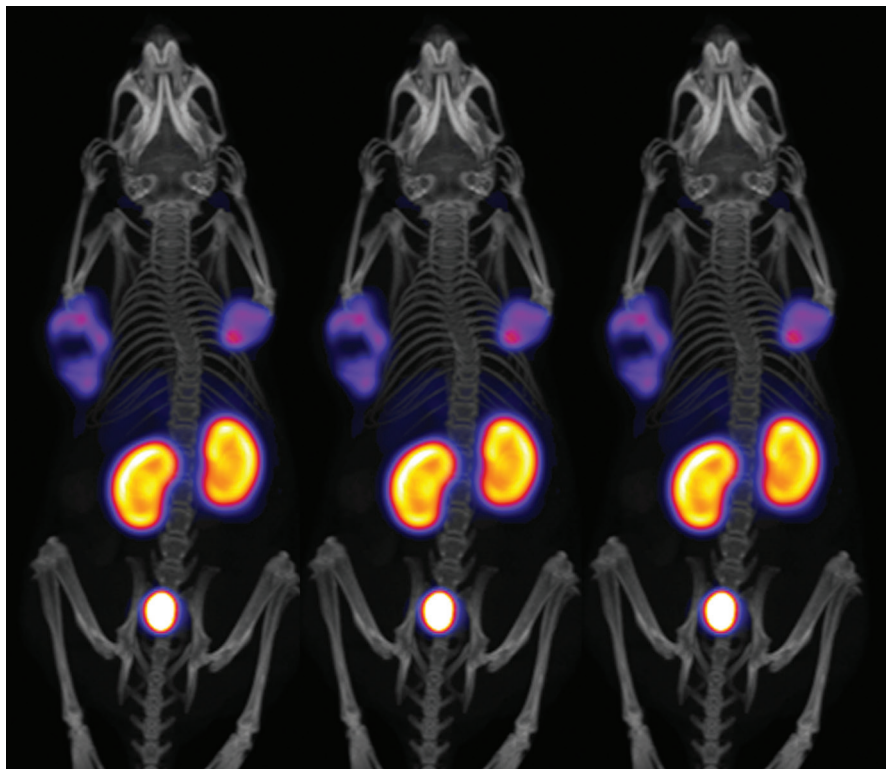


Preclinical *In Vivo* Imaging

Author:

Olivia J. Kelada, Ph.D.

PerkinElmer, Inc.
Hopkinton, MA

Radiochemistry: PET Probe Development and Imaging

Introduction

Positron Emission Tomography (PET) continues to be the primary functional imaging technique for conducting translational research studies aimed at identifying the molecular basis of human disease. This type of imaging would not be possible without the development of positron emitting radiotracers – the combination of a radionuclide with a biomolecule to target the pathway of interest. However, there are many challenges associated with radiolabeling. It is an expensive, highly-regulated and complex process that requires very specialized and inter-disciplinary knowledge. In addition, there are challenges associated with radiochemical yields, optimization of microfluidics and the chemical structures of imaging biomarkers. As a result, radiochemical yields are often low which can impact PET imaging, as generally PET scanners are not sensitive enough to detect low radiotracer activities.

With the highest sensitivity on the market, the G8 PET/CT can detect very low levels of radioactivity (Table 1), reducing the amount of activity that needs to be injected. This can alleviate the challenge of low PET probe yields and make it possible to image many biological pathways such as immune responses, where large doses of biomolecule or radioactivity could result in undesired therapeutic effects. An additional benefit is that lower amounts of probes are required to be synthesized, allowing for a more efficient use of resources, reduced procurement expenses and lower waste storage and disposal fees.

How PET Detector Sensitivity is Increased

The high PET sensitivity (~14% uniform sensitivity across the FOV) of the G8 is made possible through a) High atomic number ($Z=83$) of the BGO crystal detectors as this increases the probability of a photoelectric event at the first interaction site b) The arrangement of the four panel detector heads as they are placed close together (unlike the classic ring design) thus closer to the animal and emitted photons and c) Improved crystal cutting that reduces the gaps between adjacent crystals and increases the packing fraction. In addition, to achieve equivalent sensitivity on a

Table 1. Data demonstrates the low radiotracer activities that can be imaged on the G8 PET/CT.

Author	G4 or G8	Radiotracer	Lowest activity used (uCi)	(MBq)	Application*
Bansal et al. 2015	G4	$^{89}\text{Zr}(\text{HPO}_4)_2$, ^{89}Zr -labeled hMSCs	2.00	0.07	Novel ^{89}Zr cell labeling PET cell tracking
Gu et al. 2013	G4	^{64}Cu -minibody, ^{18}F -FDG	3.70	0.14	Performance testing of high sensitivity PET Imaging system
Yang et al. 2016	G4	$^{89}\text{Zr}(\text{HPO}_4)_2$, ^{89}Zr -labeled hMSCs	7.50	0.28	PET tracking to determine therapeutic value of human adipose tissue in AVF
Haller et al. 2015	G4	^{18}F -AzaFol, ^{18}F -fallypride	10.0	0.37	Chick and mouse comparison for PET probe uptake and stability
Meimetis et al. 2015	G4	^{89}Zr -DFO-BODIPY-trastuzumab	10.0	0.37	Multimodal imaging of antibodies with PET
Boros et al. 2016	G4	^{89}Zr -L4/L5, ^{89}Zr -DFO	10.0	0.37	Novel immunoconjugation of ^{89}Zr for PET
Peng et al. 2014	G4	^{11}C -PBR28	13.0	0.48	PET to investigate role of acetyl-coA in neurodegeneration
Lazari et al. 2014	G4	^{18}F -FLT, ^{18}F -FDG, ^{18}F -fallypride	20.0	0.74	Fully automated production of diverse ^{18}F -labeled PET tracers
Evdokimov et al. 2014	G4	^{18}F -FMT	20.0	0.74	Development of novel CNS PET probe
Faltermeier et al. 2015	G4	^{18}F -FDG	20.0	0.74	PET imaging to identify kinases driving prostate cancer metastasis
Evdokimov et al. 2015	G4	^{18}F -2-DFR, ^{18}F -DFA	20.0	0.74	Development of PET probe for liver function imaging
Kim et al. 2016	G4	^{18}F -CFA	20.0	0.74	Development of a PET probe for deoxycytidine kinase activity imaging
Hermann et al. 2013	G4	^{18}F -FDG	25.0	0.93	Evaluation of a bench-top preclinical PET scanner
Lozier et al. 2015	G4	^{18}F -FDG	25.0	0.93	Assessment of LIN28 oncogenic driver in neuroblastoma using PET
van der Meulen et al. 2015	G4	^{44}Sc -DOTANOC	27.0	1.00	Evaluation of ^{44}Sc biodistribution in vivo with PET
Boschi et al. 2016	G8	^{18}F -PSMA-11	38.0	1.40	Synthesis and preclinical evaluation of an Al ^{18}F radiofluorinated PSMA ligand
Seo et al. 2015	G4	^{64}Cu -NOTA-UCNPs	40.0	1.48	Assessment of radiolabeled upconverting nanoparticles with PET
Witney et al. 2014	G4	^{18}F -FPIA	50.0	1.85	Predicinal evaluation of a novel PET imaging agent for tumor detection
Song et al. 2016	G4	^{18}F -FMISO	50.0	1.85	PET to assess oxygen changes after high-dose radiation therapy
van Elssen et al. 2017	G8	^{64}Cu -VHH4	50.0	1.85	Imaging of human immune responses in a human xenograft model of graft vs host disease
Mueller et al. 2016	G8	^{152}Tb -DOTANOC	90.0	3.33	^{152}Tb -DOTANOC for PET radiolanthanide imaging
Farkas et al. 2016	G8	^{64}Cu - and ^{68}Ga -labeled NODAGA-folate	94.0	3.48	^{64}Cu - and ^{68}Ga -Based PET Imaging of folate receptor-positive tumors
Jang et al. 2013	G4	^{64}Cu -Sar-chTNT-3	100	3.70	PET to assess the use of chemotherapy to improve the biodistribution of antibodies
Hu et al. 2015	G4	^{18}F -FDG	100	3.70	Nanoparticle-mediated radiopharmaceutical-excited fluorescence in vivo imaging
Umbricht et al. 2017	G8	^{44}Sc -PSMA-617, ^{68}Ga -PSMA-11, ^{68}Ga -PSMA-617	135**	5.00	^{44}Sc -PSMA-617 for radiotheragnostics in comparison with ^{68}Ga -PSMA ligands
Mueller et al. 2016	G8	^{149}Tb -DOTANOC	190**	7.03	Alpha-PET with terbium-149 in mice
Domnanich et al. 2016	G8	^{44}Sc / ^{68}Ga -labeled DOTA/NODAGA peptides	270**	9.99	^{44}Sc for labeling of DOTA- and NODAGA peptides
Moon et al. 2015	G4	^{68}Ga -DOTA-IO-GUL	276**	10.2	Development of PET/MR probe for PSMA targeting

* All studies conducted in mice

** Due to tracer half-life and study design higher activities were injected prior to imaging

BGO crystal detector system, lower detector thickness is required, leading to reduced parallax errors and better spatial resolution.

Translational Research Benefits from Imaging Low Radiotracer Activities

Mice are approximately 3000 times smaller in mass and volume than humans. To obtain preclinical PET images of comparable statistical quality to clinical ones, the number of counts per voxel (which is inversely proportional to voxel volume) can either be compensated by increasing the radioactivity of injected probes or by improving the sensitivity of the scanner. However, for preclinical imaging, the typical amount of radioactivity administered is 7.4 MBq (200 μ Ci). This delivers a much higher dose per gram of tissue than in humans (typical administered activity of 370 MBq (10 mCi)). The higher absorbed dose can lead to biological effects such as stimulated cell proliferation, induced radio-resistance and elevated gene expression. Such biological effects may interfere and bias the results in pharmaceutical and genetic studies, leading to discrepancies in translational research between mouse models and clinical applications. This may be even more important for longitudinal studies that require serial imaging. Also, for many applications, signal uptake can be limited by the number of binding sites available, thus when higher activity of radiotracer (and in turn higher amounts of targeting agent) are injected, this may block the binding receptors and prevent repeated imaging. The amount of injected probe should be below levels that perturb the in vivo system and this may limit the amount of radioactivity that can be injected into the animal. Thus,

a preclinical PET scanner with high sensitivity lowers the activities of radioisotopes that are needed to obtain high quality PET images. In turn, this results in less exposure to the animal and the investigator. For biodistribution studies on the G8, the typical administered radioactivity amount should be less than 1.5 MBq (40 μ Ci), which is five times lower than established procedures and is potentially more physiologically relevant¹.

Novel Radiotracer Development Using the G8 PET/CT

Figure 1 shows images of newly developed tracers using the G8 PET/CT.

Panel a ^{64}Cu -rf42 (an albumin binding NODAGA-folate) showed much better resolution than ^{68}Ga -rf42 of tumor xenografts and kidneys in CD1 nude mice with KB tumor xenografts and SCID mouse with a LNCaP tumor xenografts. The image shows a 10 min static scan of 4.3 MBq (116.2 μ Ci) and standard CT. The data suggests that ^{64}Cu -rf42 is a more favorable PET agent due to improved tumor-to-kidney ratios and the more physiologically suitable half-life of ^{64}Cu as it matches the circulation time of the albumin binding NODAGA-folate and allows for serial PET imaging at later time points which is not possible when using ^{68}Ga ².

Panel b ^{18}F -AzaFol (^{18}F -labeled folate-based radiotracer) shows specificity in both the tumor and kidneys in CD1 nude mice with KB tumor xenografts. The image shows a 10 min static scan of 2.1 MBq (56.8 μ Ci) and standard CT. Results suggested that ^{18}F -AzaFol is an appropriate tool for imaging FR-positive tumors³.

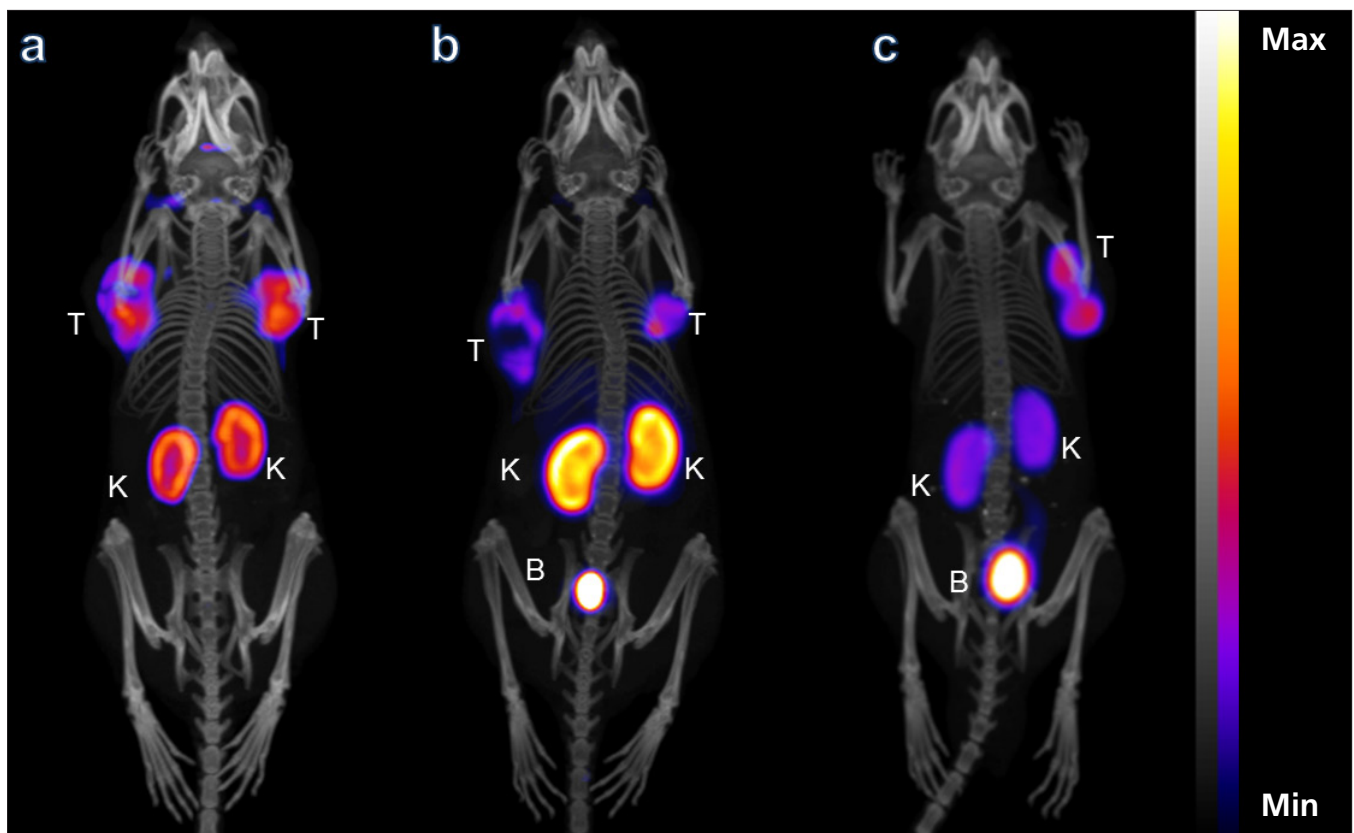


Figure 1. ^{64}Cu -NODAGA-folate static PET scan in CD1 nude mice with cervical cancer xenografts (b) ^{18}F -AzaFol static PET in CD1 nude mice with cervical cancer xenografts (c) ^{44}Sc -labeled PSMA-617 static PET/CT scan of SCID mouse with LNCaP prostate cancer xenograft T = Tumor, K = Kidney and B = Bladder All PET/CT images courtesy of Dr. Cristina Müller Center for Radiopharmaceutical Sciences (ETH/PSI), Zurich, Switzerland and used with permission.



PET Probe Development and Imaging on the G8 PET/CT: Customer Testimonial

Dr. Cristina Müller Center for Radiopharmaceutical Sciences (ETH/PSI), Zurich, Switzerland

"The G8 PET/CT was installed at the Center for Radiopharmaceutical Sciences ETH-PSI in 2015. It has proven to be particularly effective for the evaluation of novel in-house produced radiotracers, that are initially only available in small quantities. The fact that the scanner is small and mobile has allowed us to transport it to other facilities including the Center for Proton Therapy located at the PSI where the G8 scanner was used for imaging in vivo ^{11}C production after proton irradiation of tumor xenografts in mice."

Panel c ^{44}Sc labeled PSMA-617 shows specificity and sensitivity in SCID mouse with LNCaP tumor xenografts. The image shows a 10 min static scan of 0.8 MBq and standard CT. Due to the desirable nearly four-fold longer half-life of ^{44}Sc compared to ^{68}Ga , the feasibility of centralized production of ^{44}Sc and transport to satellite PET centers, ^{44}Sc -PSMA-617 is appealing for clinical PET imaging of prostate cancer patients⁴.

Summary

With the highest sensitivity on the market, the preclinical G8 PET/CT can detect very low levels of radioactivity. This is ideal for PET probe development, where yields are often low, and studies investigating various applications where large doses of carrier or radioactivity could elicit unexpected biological effects or block the binding sites and prevent repeated imaging. The amount of injected probe should be below levels that perturb the in vivo system and this may limit the amount of radioactivity that can be injected into the animal. The G8 PET/CT facilitates imaging with more translation doses. This, in combination with its small footprint, makes it the ideal tool for the advancement of novel PET radiotracer development and imaging.

References

1. Gu Z, et al. NEMA NU-4 performance evaluation of PETbox4, a high sensitivity dedicated PET preclinical tomograph. *Phys Med Biol.* 2013; 58:3791-814.
2. Farkas R, et al. ^{64}Cu - and ^{68}Ga -Based PET Imaging of Folate Receptor-Positive Tumors: Development and Evaluation of an Albumin-Binding NODAGA-Folate. *Mol. Pharmaceutics* 2016, 13,1979-87.
3. Betzel T, et al. Radiosynthesis and Preclinical Evaluation of 3'-Aza-2'-[^{18}F]fluorofolic Acid: A Novel PET Radiotracer for Folate Receptor Targeting. *Bioconjug Chem* 2013, 24, 205-14.
4. Umbricht CA, et al. ^{44}Sc -PSMA-617 for radiotheragnostics in tandem with ^{177}Lu -PSMA-617 - preclinical investigations in comparison with ^{68}Ga -PSMA-11 and ^{68}Ga -PSMA-617. *EJNMMI Research* 2017, 7:9.