

DIPG Model

- Diffuse intrinsic pontine glioma (DIPG) is an aggressive pediatric tumor of the pons, responsible for critical autonomic and motor functions. It accounts for 80% of childhood brainstem tumors and typically affects children age 5–10, with approximately 300 U.S. cases per year. The tumor is highly infiltrative and unresectable; radiation offers only transient benefit, and median survival remains 9–12 months.
- Subcutaneous (s.c.) implantation shows poor tumor growth in both NCG and Athymic Nude mice, limiting utility for standard s.c. in vivo studies.
- Orthotopic implantation into the pons results in tumor growth, consistent with DIPG biology.
- Pons-resident tumors maintain disease-relevant behavior, supporting use for translational studies.

DIPG *in vivo* Growth

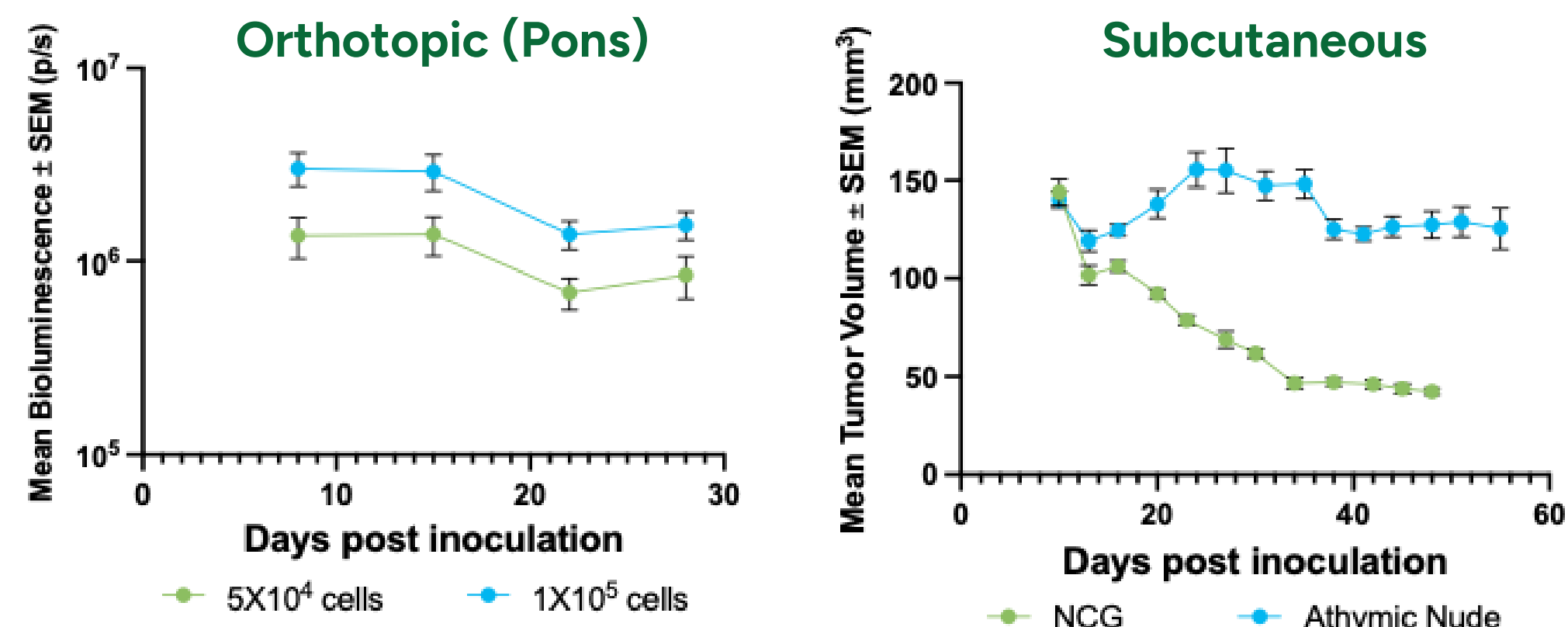


Figure 1: Comparison of SF8628 tumor growth following pons versus subcutaneous inoculation. A localized luminescent signal was observed in the pons region of athymic nude mice receiving intracranial inoculations of 5×10^4 or 1×10^5 cells. In contrast, subcutaneous inoculation of 1×10^7 cells (1:1 Matrigel:Serum-Free Media) in NCG mice resulted in tumor regression to volumes below 50 mm^3 over time. Subcutaneous implantation of 4×10^6 cells (1:1 Matrigel:Serum-Free Media) in athymic nude mice produced tumors that stabilized at approximately 120 mm^3 .

Acknowledgements



The MRI Core (RRID:SCR_021198) services and instruments used in this project were funded, in part by the Pennsylvania State University College of Medicine via the Office of the Vice Dean of Research and Graduate Students and the Pennsylvania Department of Health using Tobacco Settlement Funds (CURE). The content is solely the responsibility of the authors and does not necessarily represent the official views of the university or College of Medicine. The Pennsylvania Department of Health specifically disclaims responsibility for any analyses, interpretations, or conclusions.



Contact Information

Cheryl Davis, MLAS, rLATg

Vice President US In Vivo



Reaction Biology Hershey
1214 Research Boulevard
Suite 1050
Hummelstown, PA 17036



+1 717-910-1437



Cheryl.Davis@reactionbiology.com

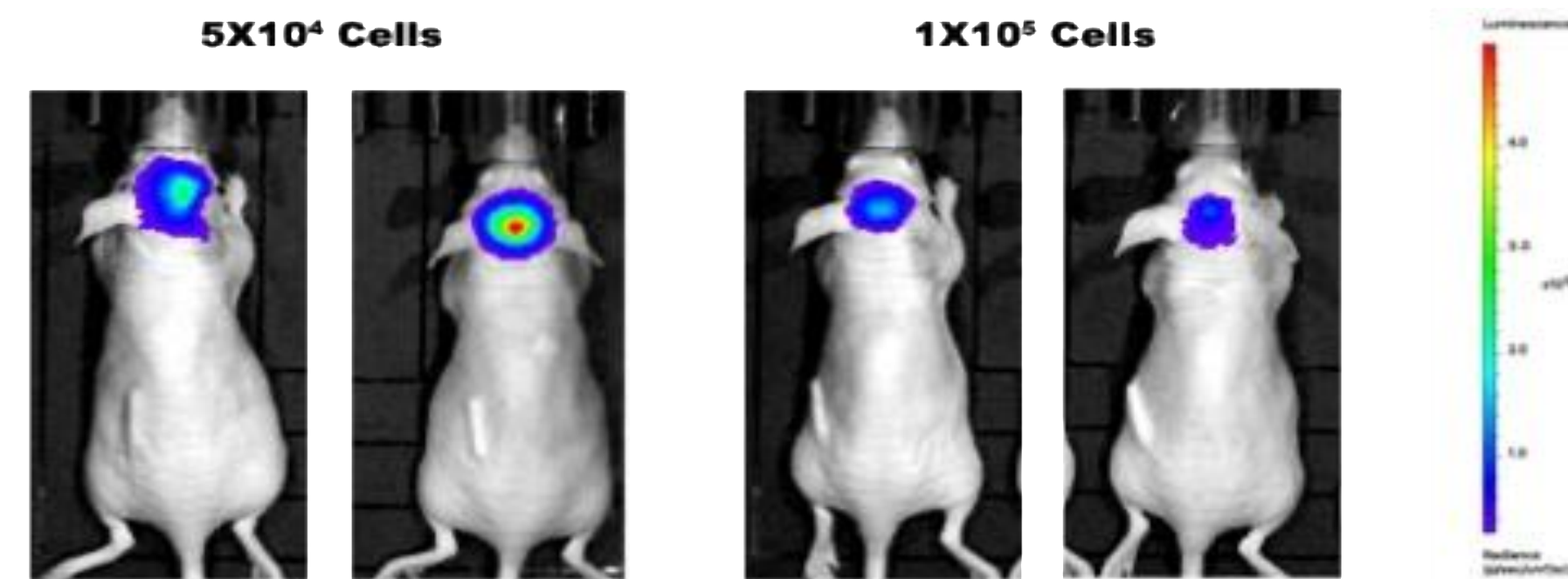


www.reactionbiology.com

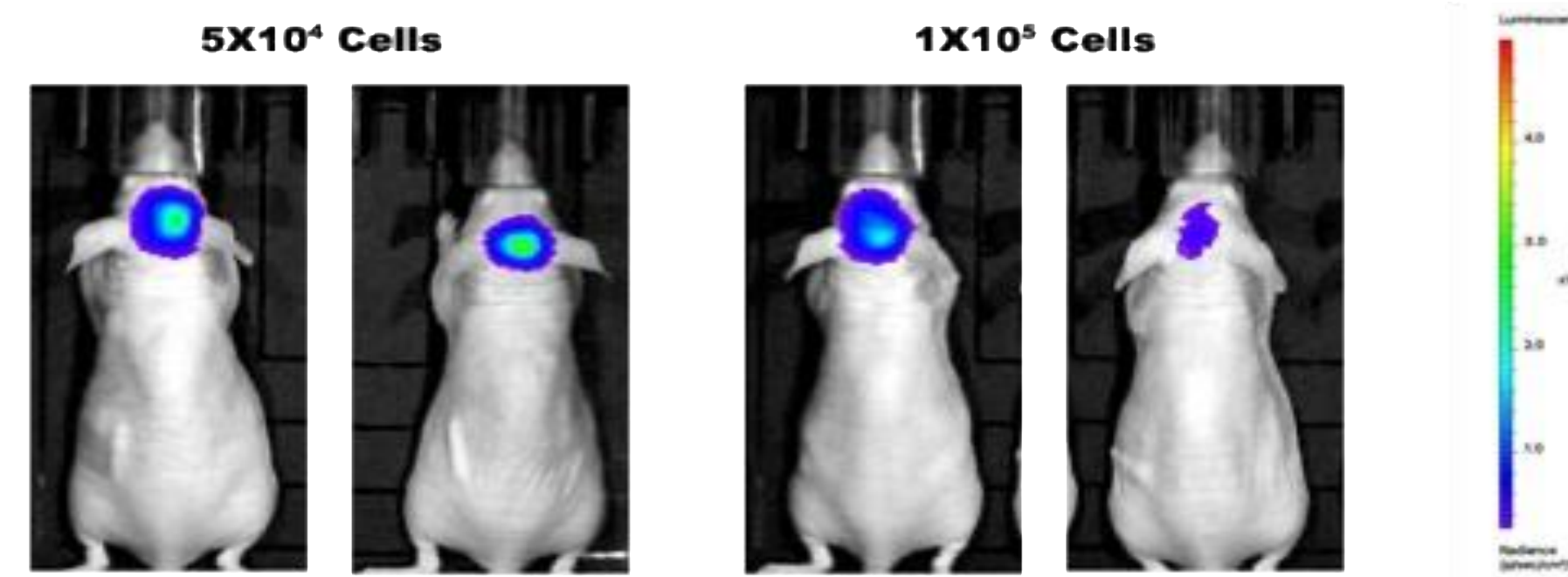
DIPG In-Life Imaging

Bioluminescent Imaging (BLI)

8 Days Post Inoculation



28 Days Post Inoculation



Magnetic Resonance Imaging (MRI)

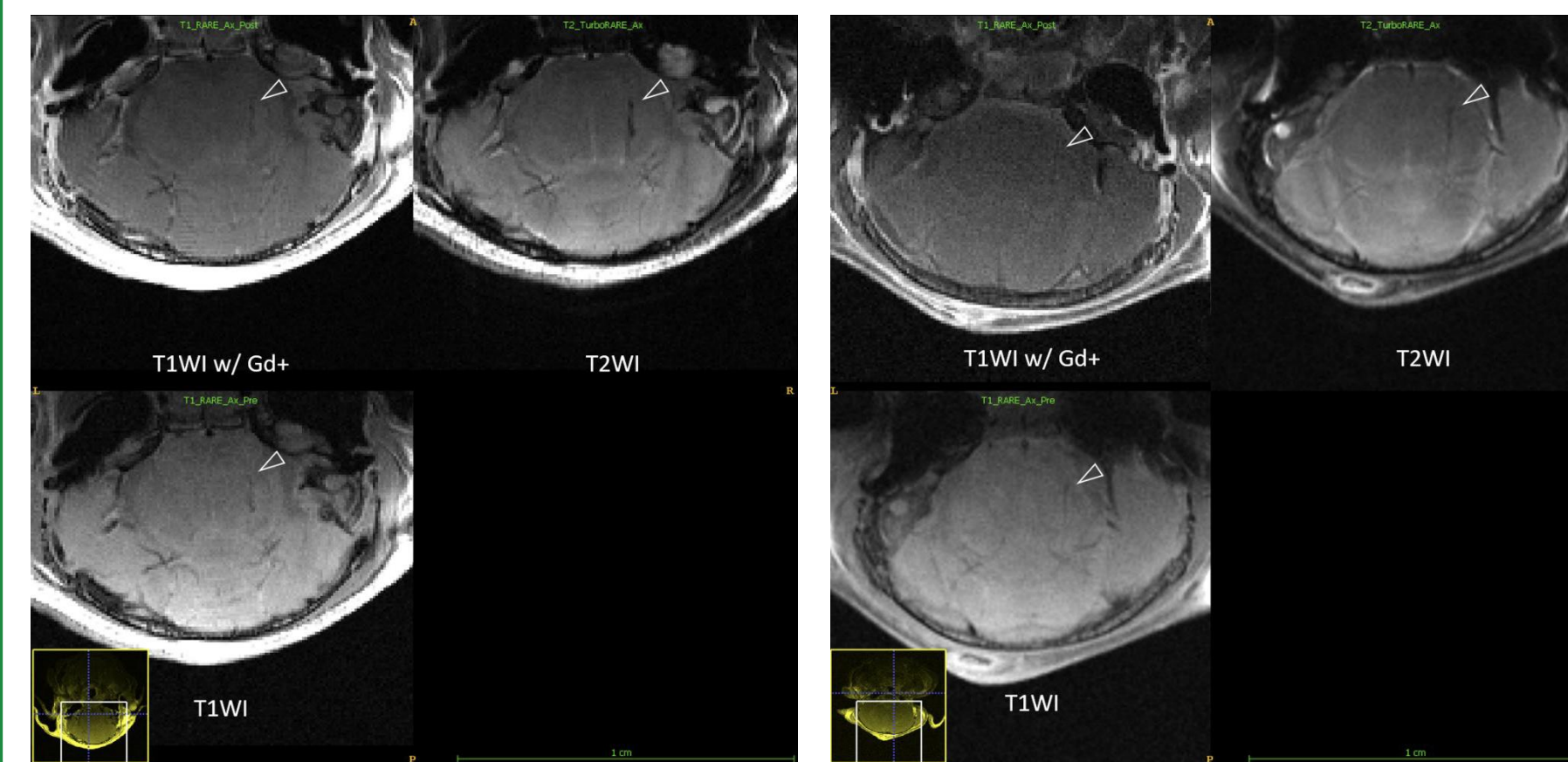


Figure 2: Evaluation of DIPG Tumor Growth in the Pons. Luminescent signal from pons-inoculated DIPG tumor cells was detectable at Day 8 and remained present at Day 28, with a modest decline possibly due to reduced luciferin delivery, hypoxia, and photon attenuation. Thirty days post inoculation, MRI showed a 5.4 mm inoculation tract extending from the skull surface to the right pons (short T2-weighted signal ~1.2 mm from the midline). A small subcutaneous chronic hemorrhage near the injection site was visible in both animals ($2.5 \times 0.2 \text{ mm}$ and $1.2 \times 0.4 \text{ mm}$).

Meningioma Model

- Meningiomas are the most common primary central nervous system tumors, constituting nearly 40% of intracranial tumors, and while generally benign, some exhibit invasive growth and recurrence requiring repeated treatment.
- Subcutaneous implantation models often show poor or unreliable tumor development, limiting their usefulness for translational studies, whereas orthotopic intracranial or subdural implantation produces tumors that closely mimic human meningiomas.
- Subcutaneous (s.c.) implantation shows poor tumor growth in C.B-17 scid mice, limiting utility for standard s.c. in vivo studies.
- Orthotopic implantation into the meningioma results in tumor growth, consistent with meningioma biology.
- Subdural-resident tumors maintain disease-relevant behavior, supporting use for translational studies.

Meningioma *In Vivo* Growth

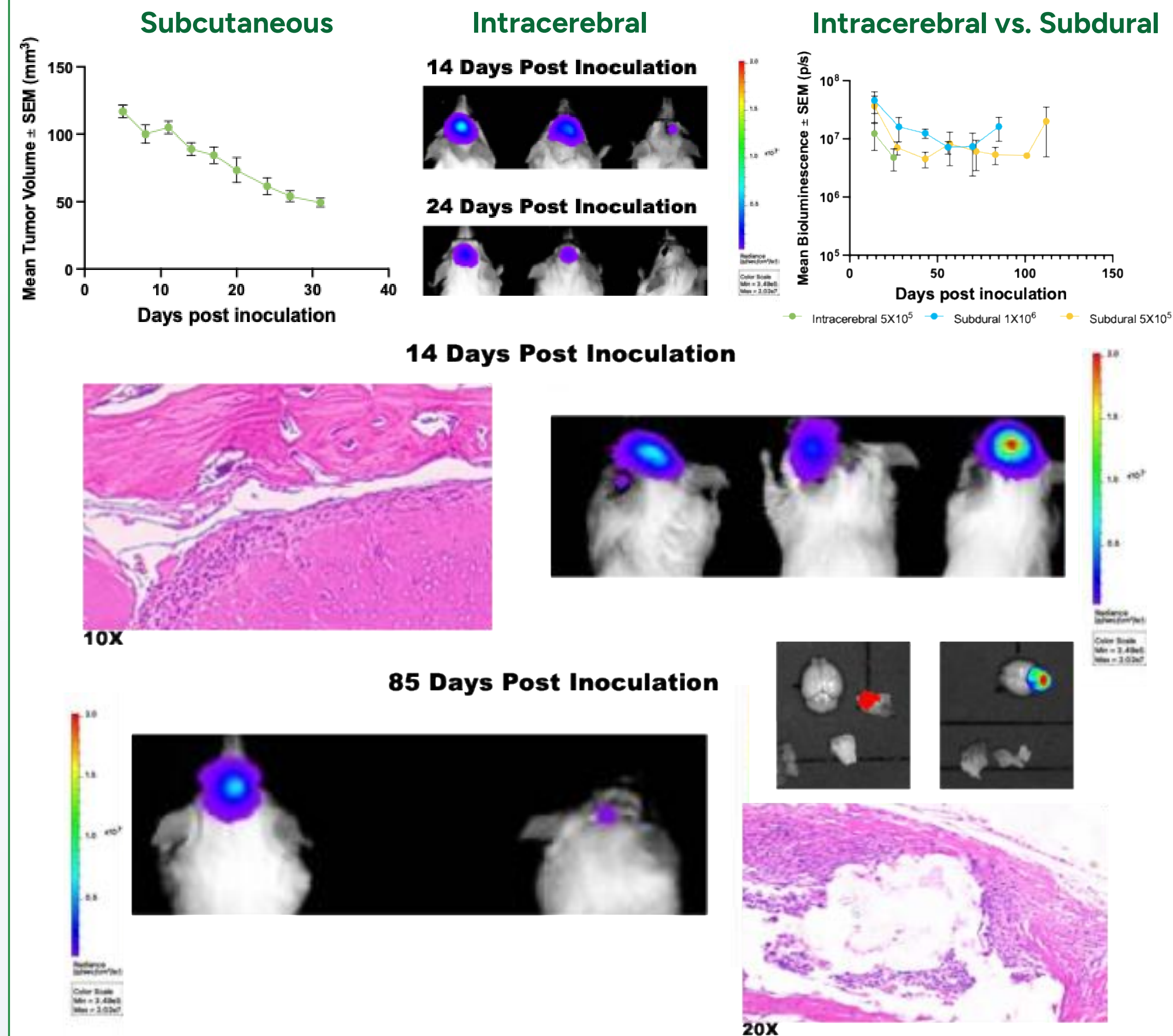


Figure 3: Comparison of Ben Men-1-luc tumor growth in C.B-17 SCID mice following subcutaneous, intracerebral, or subdural implantation. Subcutaneous inoculation of 1×10^7 Ben Men-1-luc cells (1:1 Matrigel:Serum-Free Media) resulted in declining tumor size over time, with an average tumor volume of 49.34 mm^3 at Day 31. Intracerebral implantation of 5×10^5 cells (1:1 Matrigel:HBSS) produced a luminescent signal that decreased by Day 24. In contrast, subdural implantation of 5×10^5 or 1×10^6 cells generated stable bioluminescent signal from Days 27 and 21 onward, respectively, with evidence of progressive growth by Days 101 and 85. *Ex vivo* luminescent and H&E imaging at termination confirmed tumors localized either to the brain surface or inner skull.

References

- Diffuse Intrinsic Pontine Glioma – UpToDate (Marcus et al., 2025).
- Khan M, et al., Modeling Meningiomas, Neurosurgery Clinics, 2023; 34, 479-492