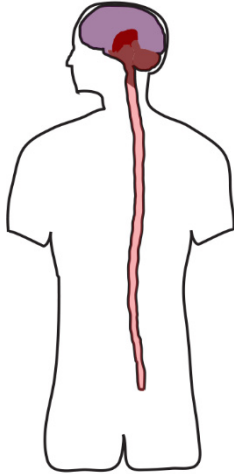


Optical control of neural ablation in zebrafish as a model for secondary injury mechanisms

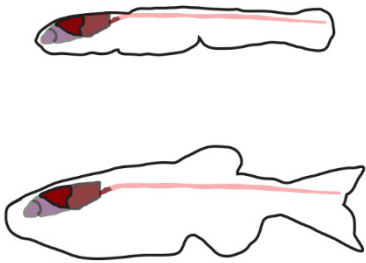
Karen Mruk, Patrick A. Piza, and James K. Chen
Departments of Chemical and Systems Biology,
Developmental Biology, and Chemical Engineering

Stanford, University, Stanford, CA

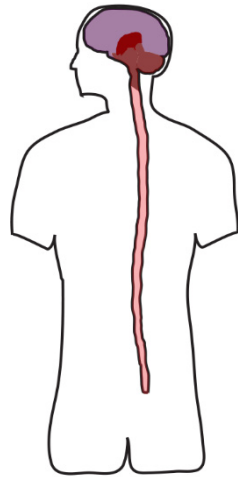
Behavioral consequences of neuronal loss



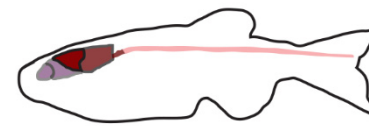
Complex Vertebrates
Myelinated Axons
Neurotransmitters



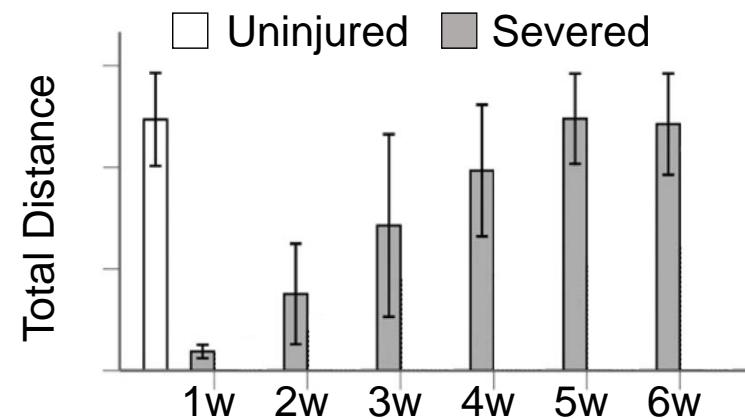
Behavioral consequences of neuronal loss



Complex Vertebrates
Myelinated Axons
Neurotransmitters



Permanent
Dysfunction



Adapted from: Yu, Y and Schachner, M. *Eur J Neurosci.* 2013

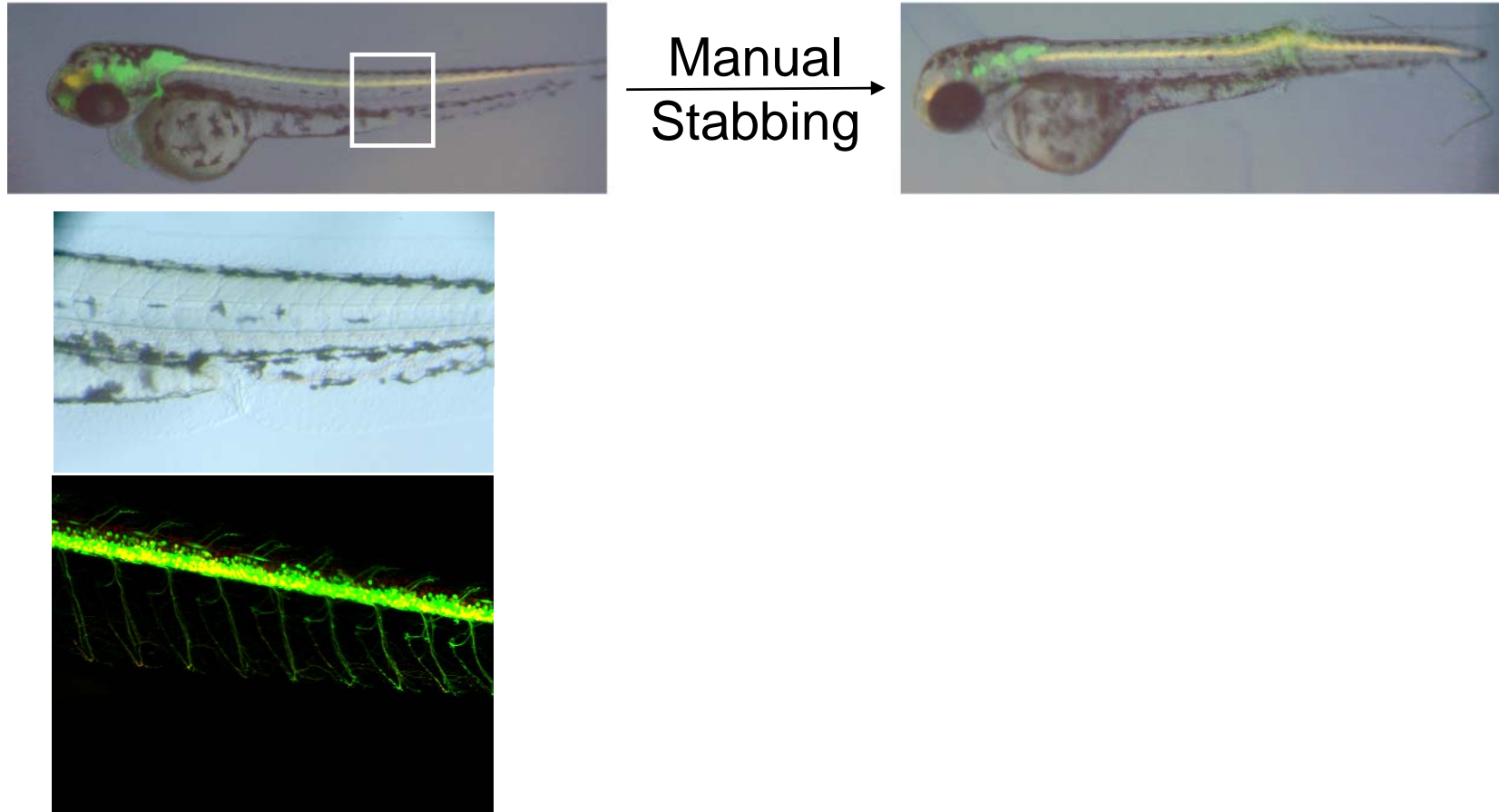
Current models of spinal injury in zebrafish are limited



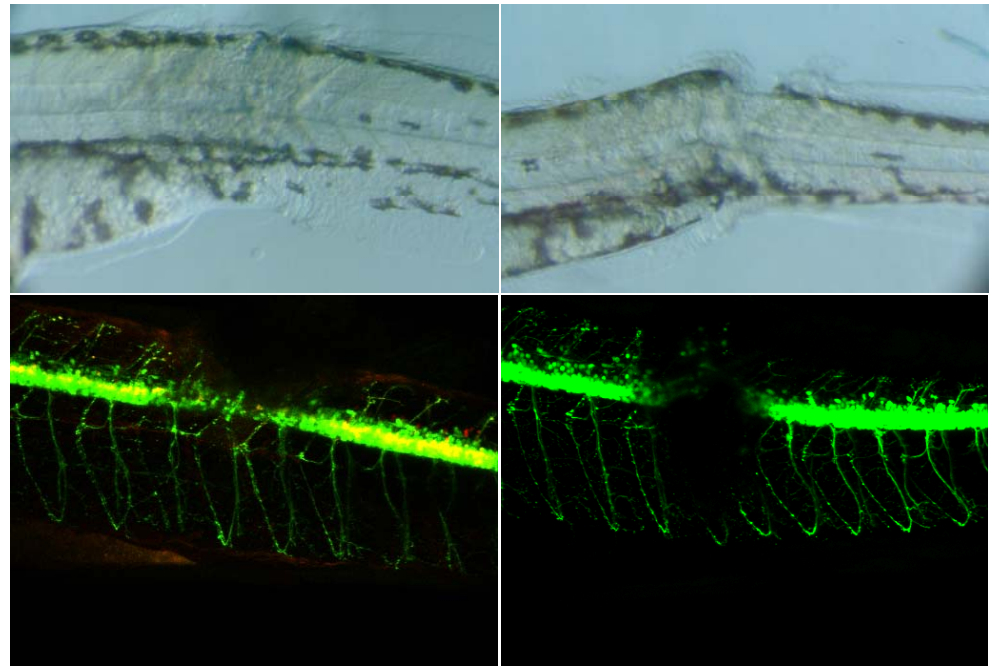
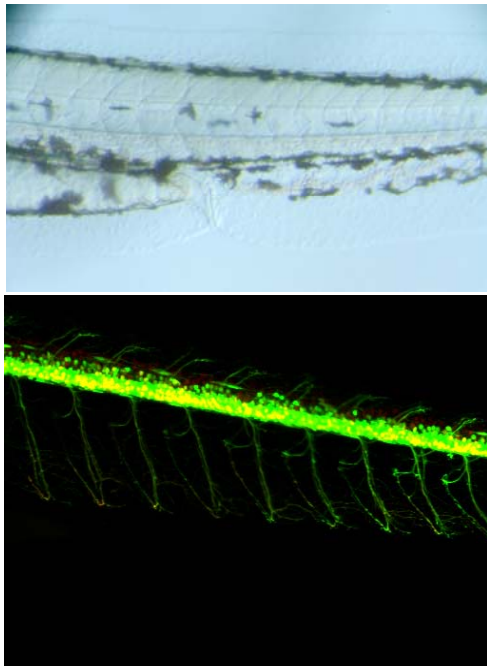
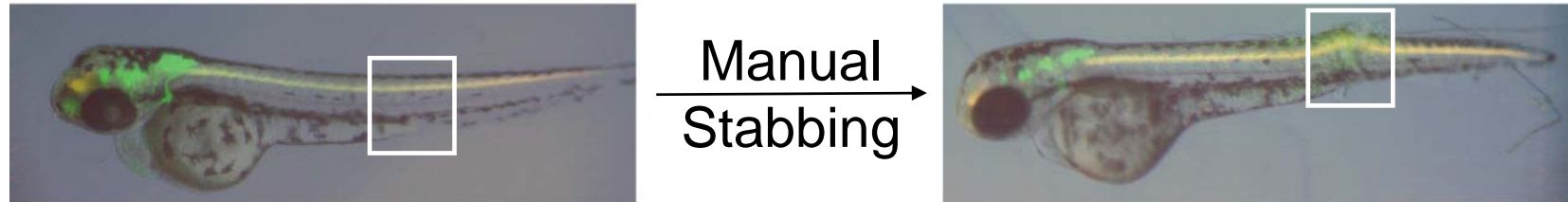
Manual
Stabbing →



Current models of spinal injury in zebrafish are limited

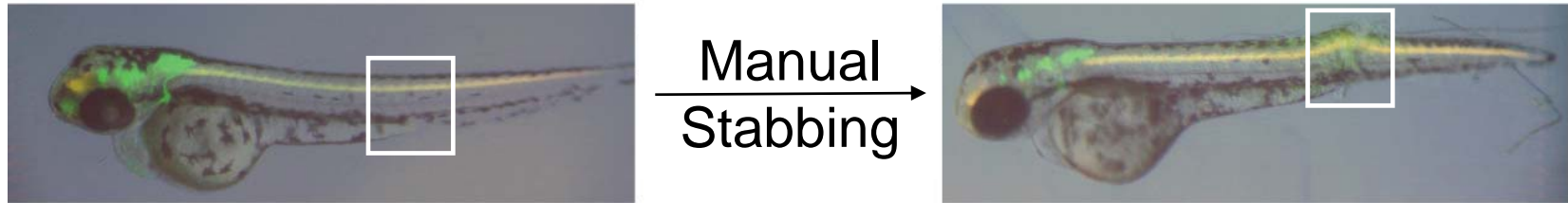


Current models of spinal injury in zebrafish are limited



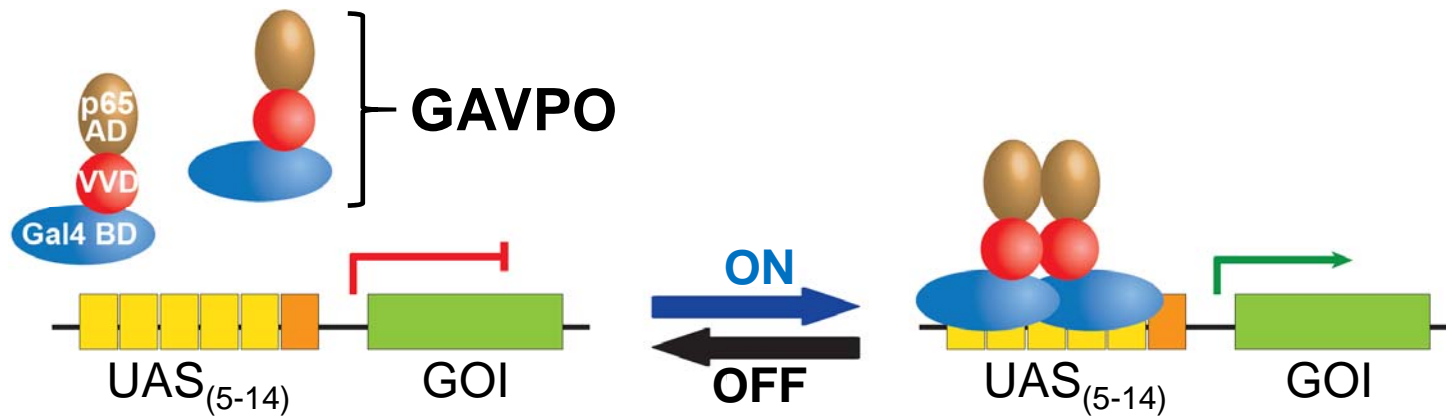
Less reproducible
Extraneous damage
Higher mortality

Current models of spinal injury in zebrafish are limited



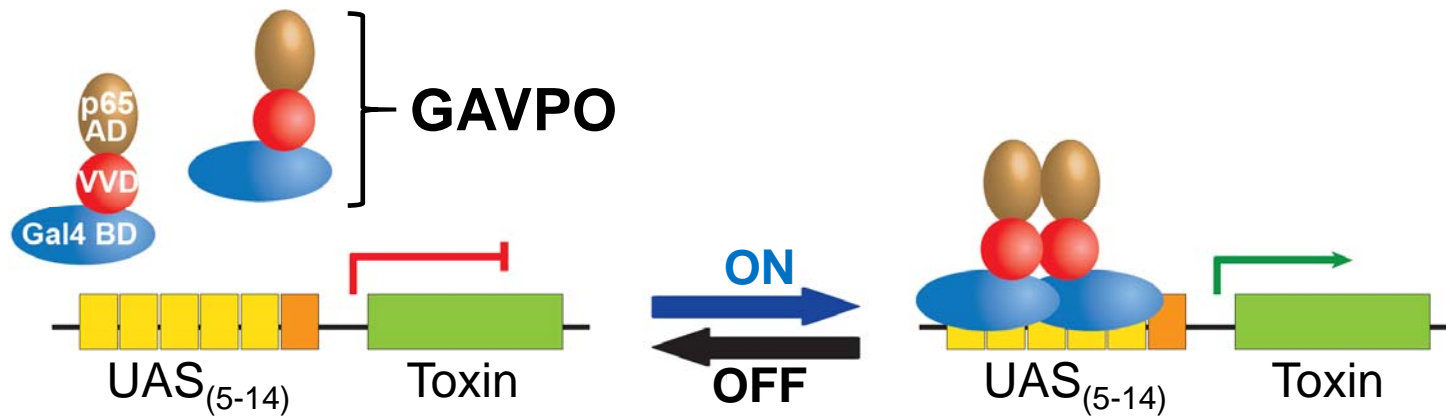
Can we use an optogenetic approach to reproducibly kill neurons as a model for spinal cord injury and degeneration?

Using light to spatiotemporally ablate CNS cells

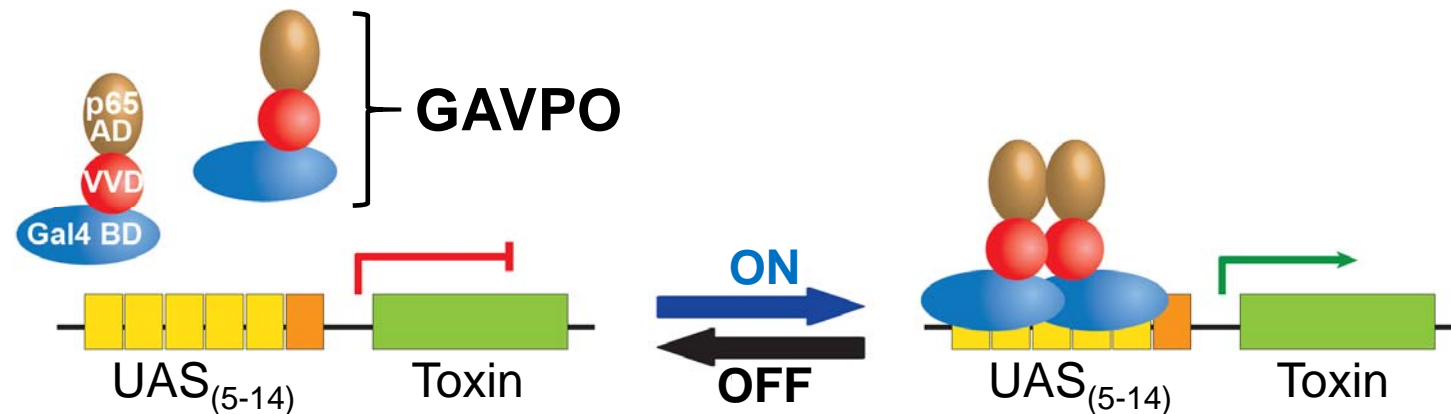


Wang et al. *Nature Methods* 2012

Using light to spatiotemporally ablate CNS cells

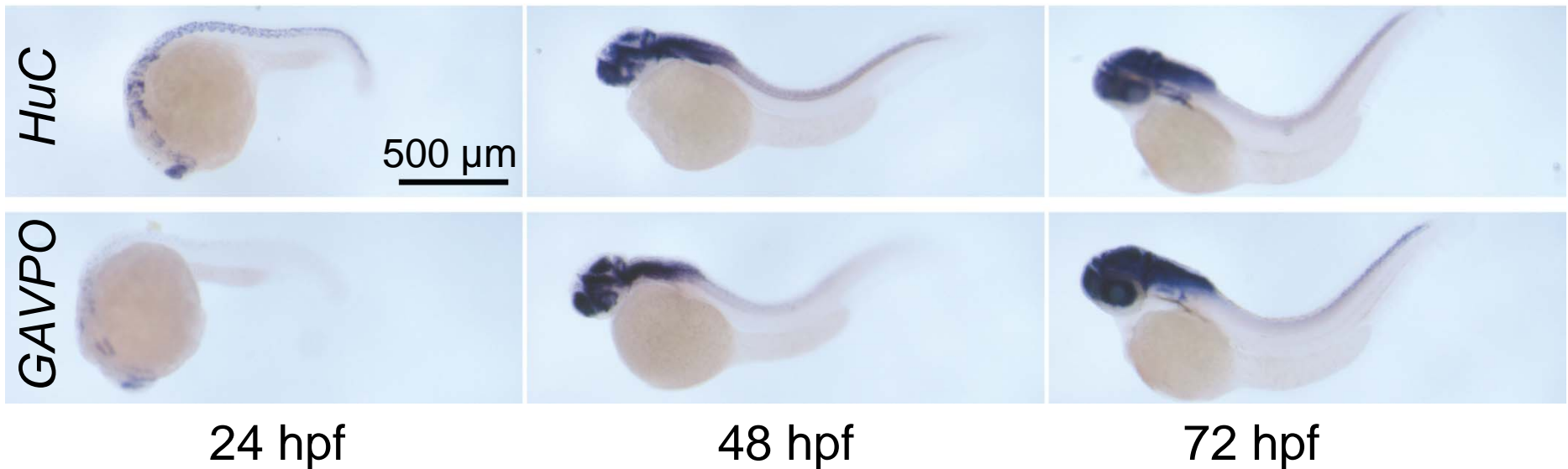
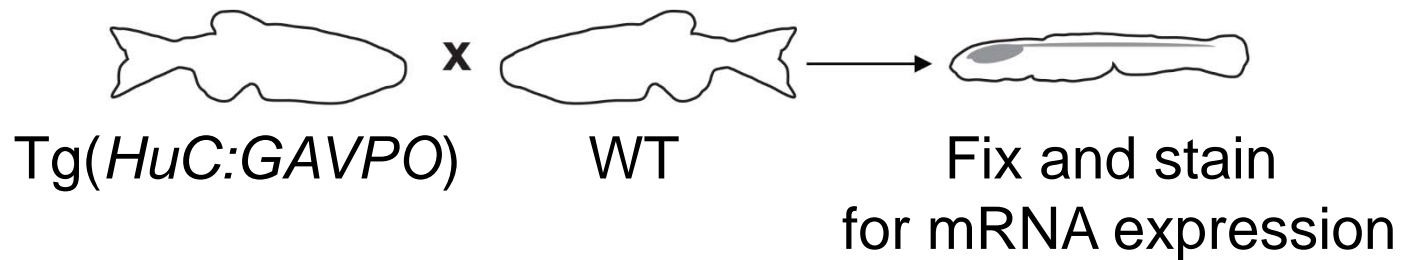


Using light to spatiotemporally ablate CNS cells



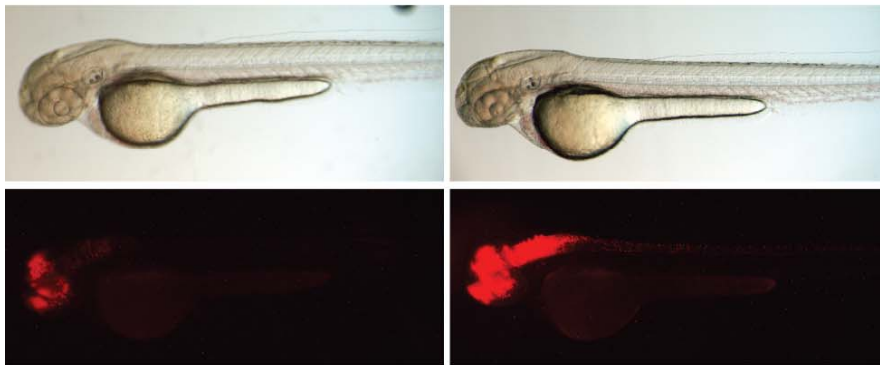
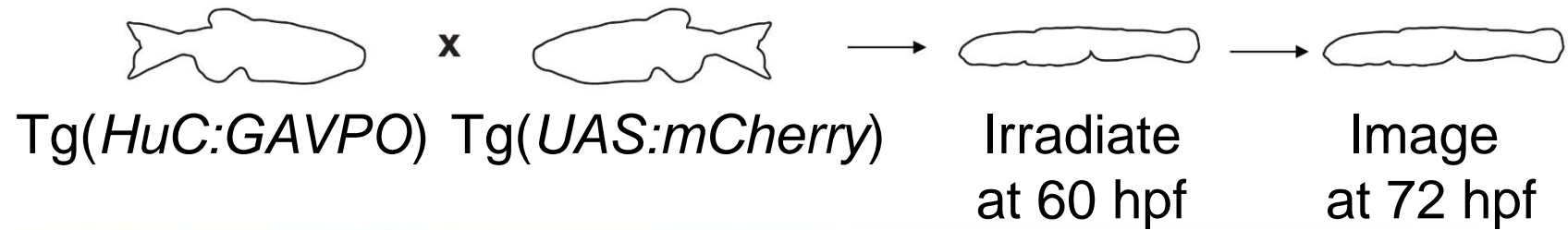
- Does GAVPO drive transcription in zebrafish?
- Can we model neural death with toxins?
- Can we induce neural death using the GAVPO system?

Generating and characterizing transgenic GAVPO lines



n = 20-25 larvae for each condition

GAVPO-induced expression increases with irradiation duration and intensity



20 min

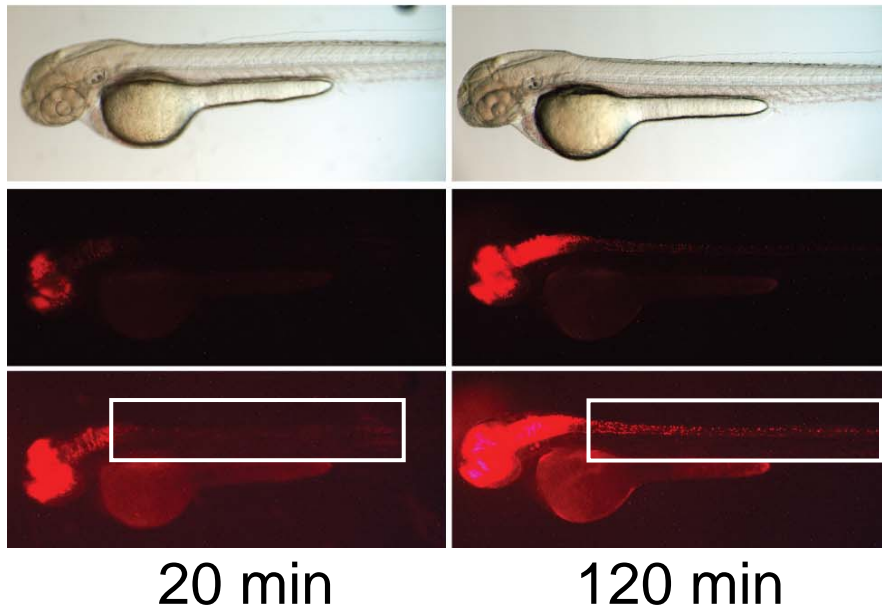
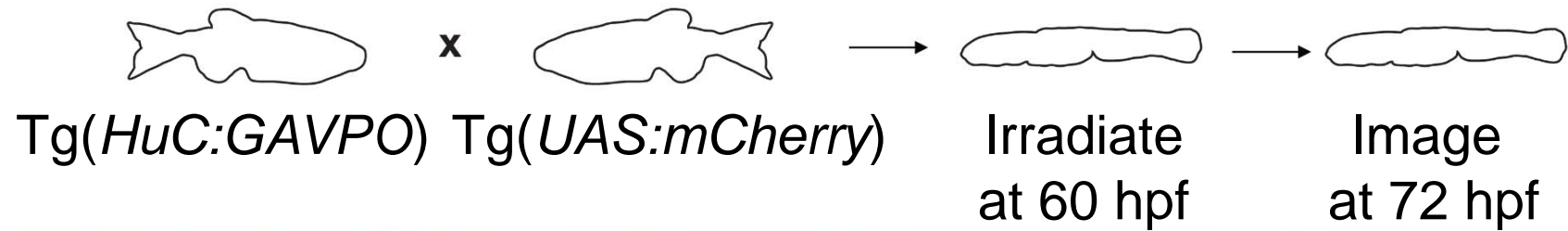
120 min

Global Irradiation

3 mW/cm²

n = 25-30 larvae for each condition

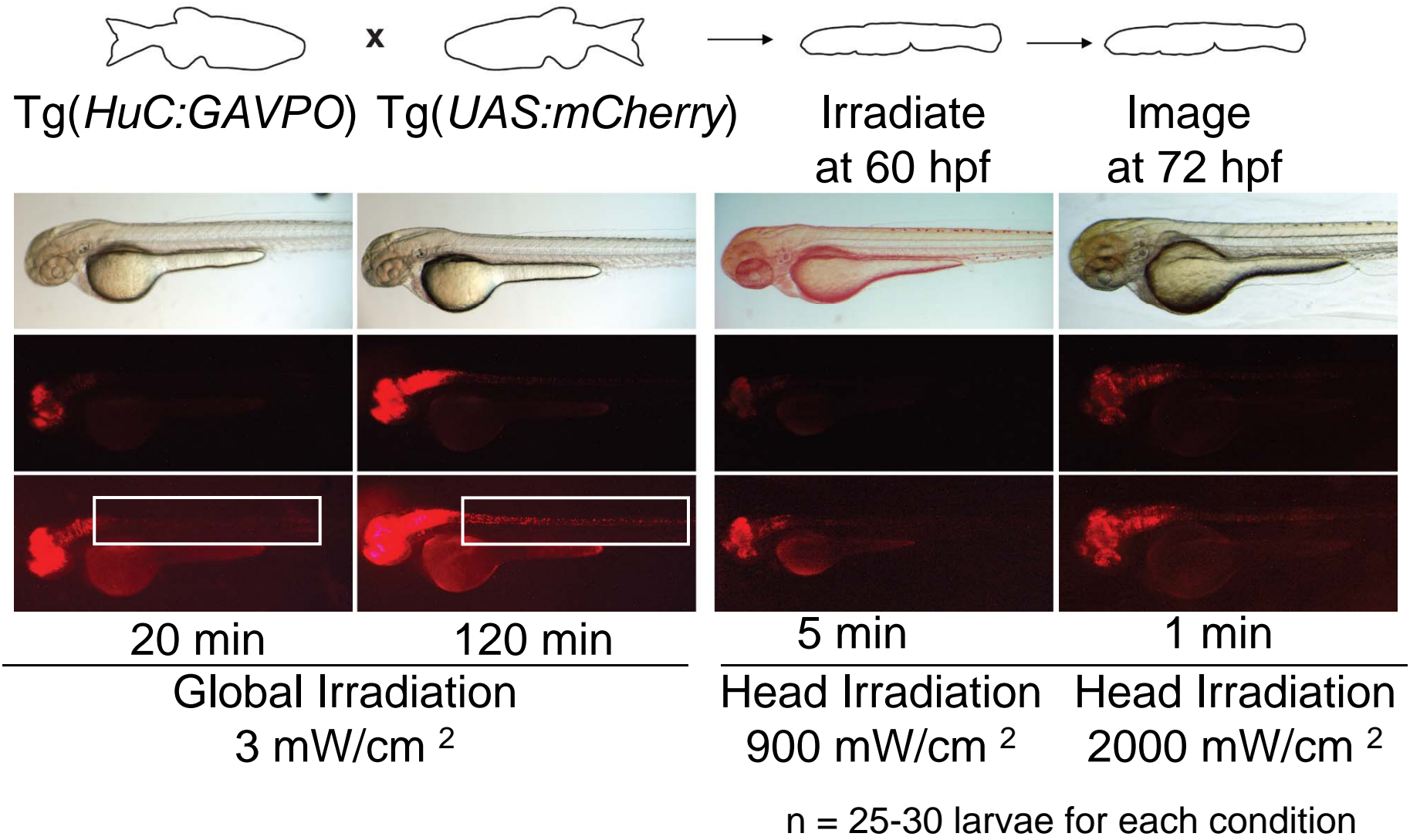
GAVPO-induced expression increases with irradiation duration and intensity



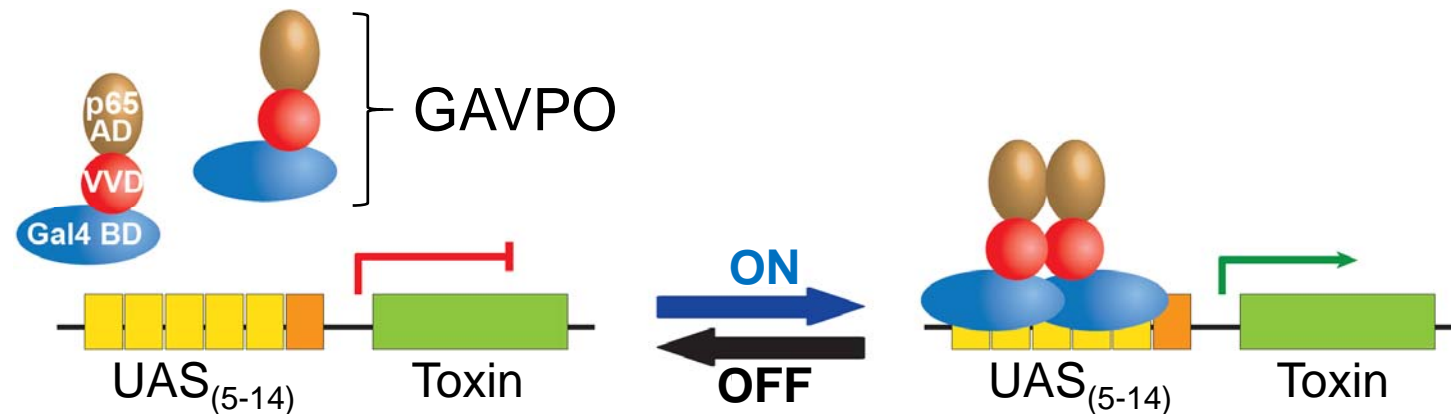
Global Irradiation
3 mW/cm²

n = 25-30 larvae for each condition

GAVPO-induced expression increases with irradiation duration and intensity

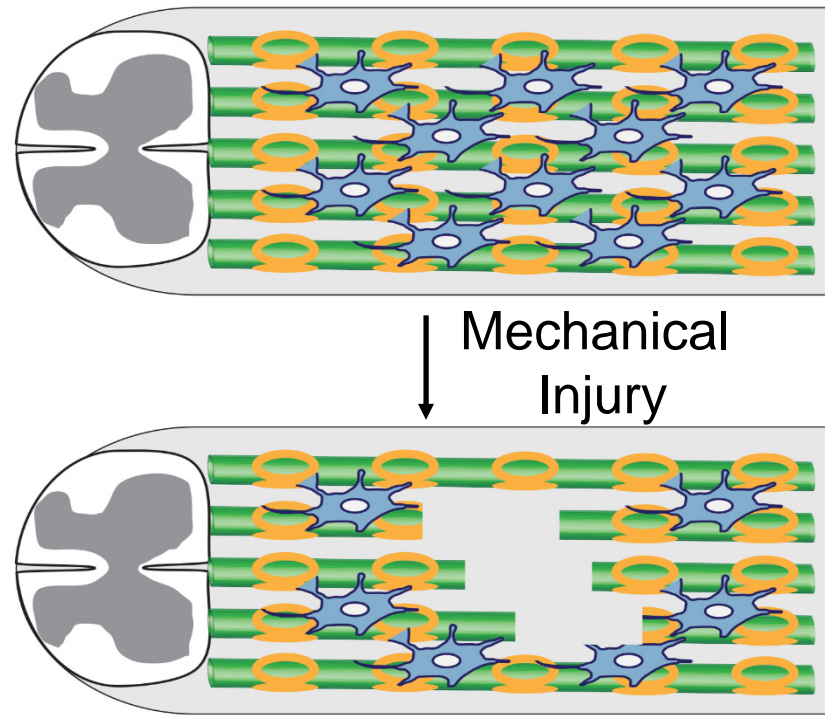


Using light to spatiotemporally ablate CNS cells

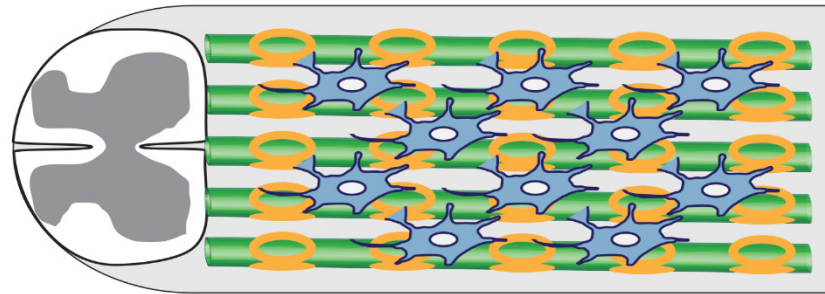


- ✓ Does GAVPO drive transcription in zebrafish?
- ❑ Can we model neural death with toxins?
- ❑ Can we induce neural death using the GAVPO system?

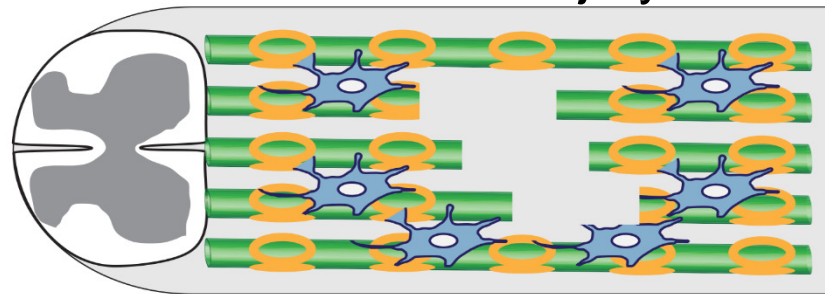
Modeling secondary injury via toxin-mediated ablation



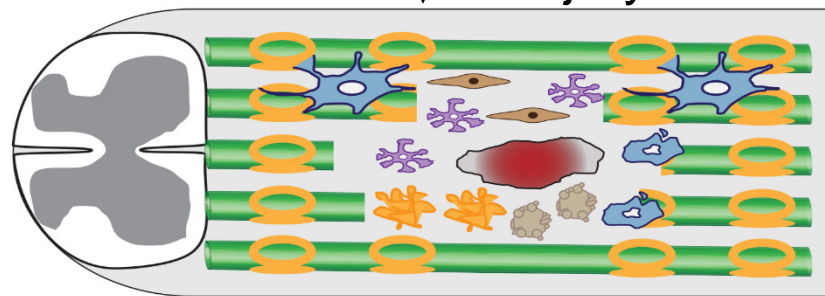
Modeling secondary injury via toxin-mediated ablation



Mechanical Injury

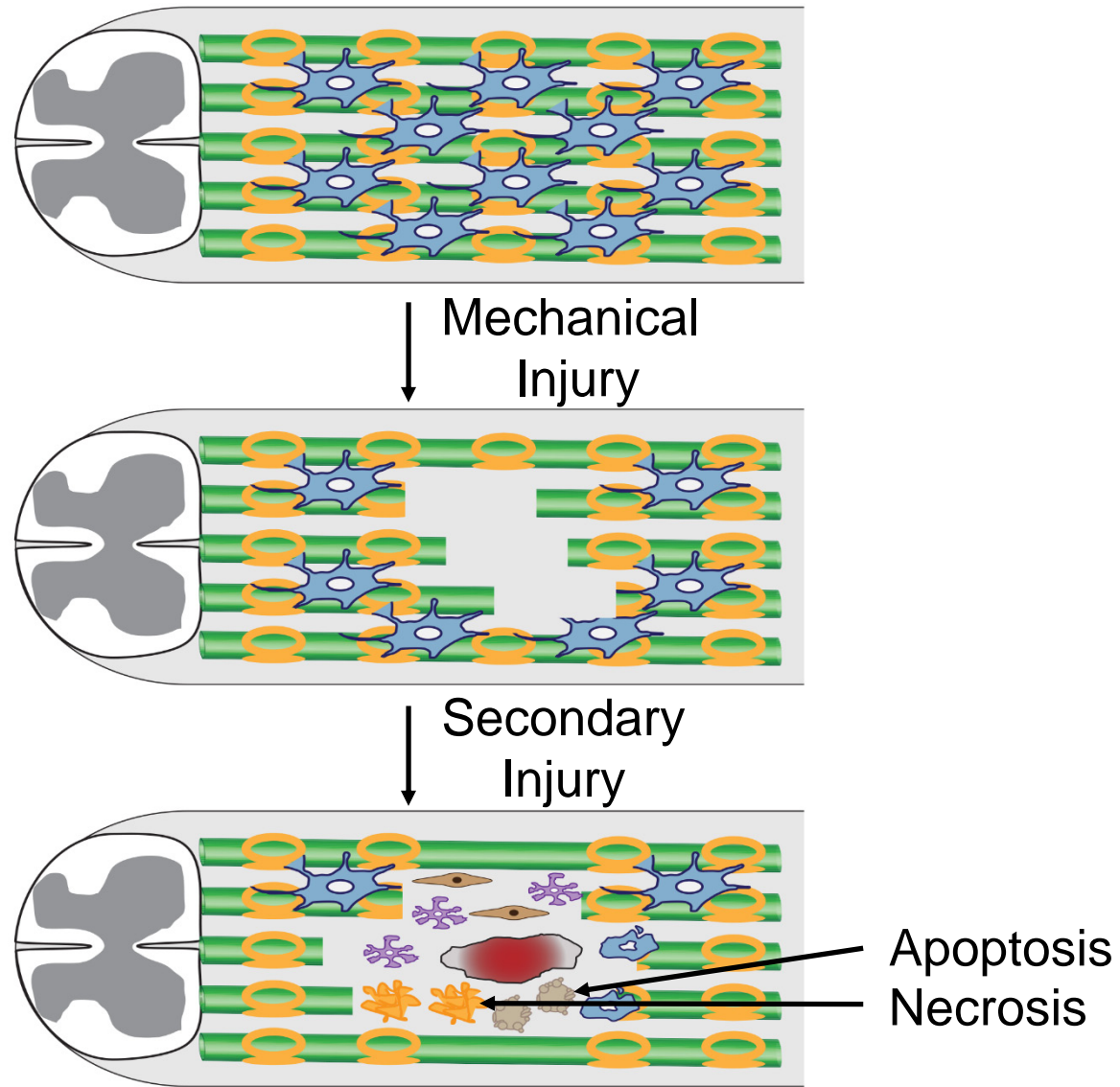


Secondary Injury

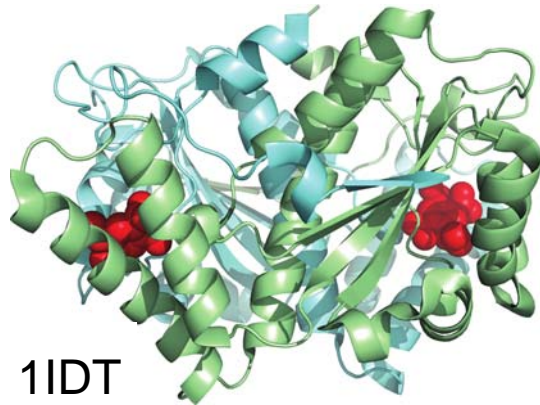


Cell infiltration
ROS generation
Cell death

Modeling secondary injury via toxin-mediated ablation



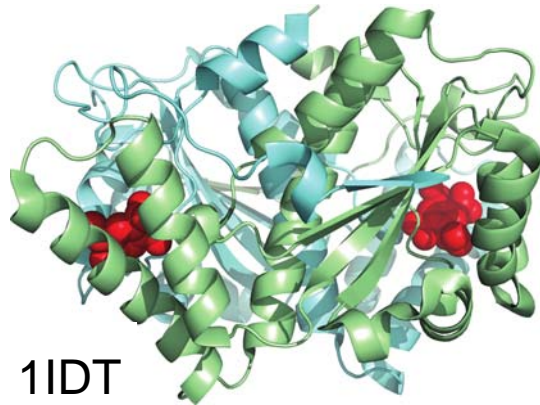
Modeling neuronal loss via development of toxin-mediated ablation



PDT: 1IDT

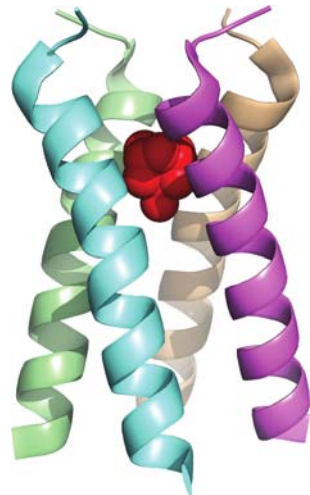
- Nitroreductase
 - Enzyme converts a prodrug (metronidazole) into a cytotoxic compound
 - Gold standard for zebrafish
 - **Disadvantage – slow acting**

Modeling neuronal loss via development of toxin-mediated ablation



PDT: 1IDT

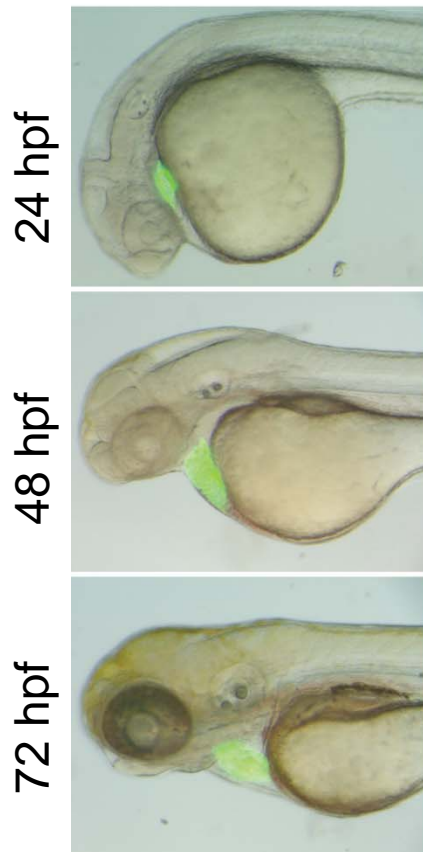
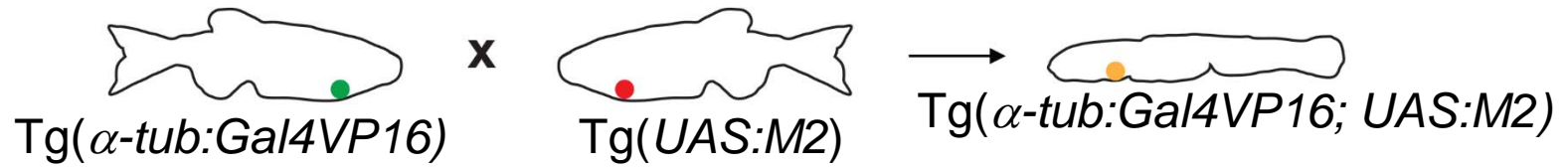
- Nitroreductase
 - Enzyme converts a prodrug (metronidazole) into a cytotoxic compound
 - Gold standard for zebrafish
 - **Disadvantage – slow acting**



PDT: 2KQT

- M2
 - Ion channel from flu
 - Can be blocked with FDA approved drug (rimantadine)
 - Never used in zebrafish
 - **Advantage – fast acting**

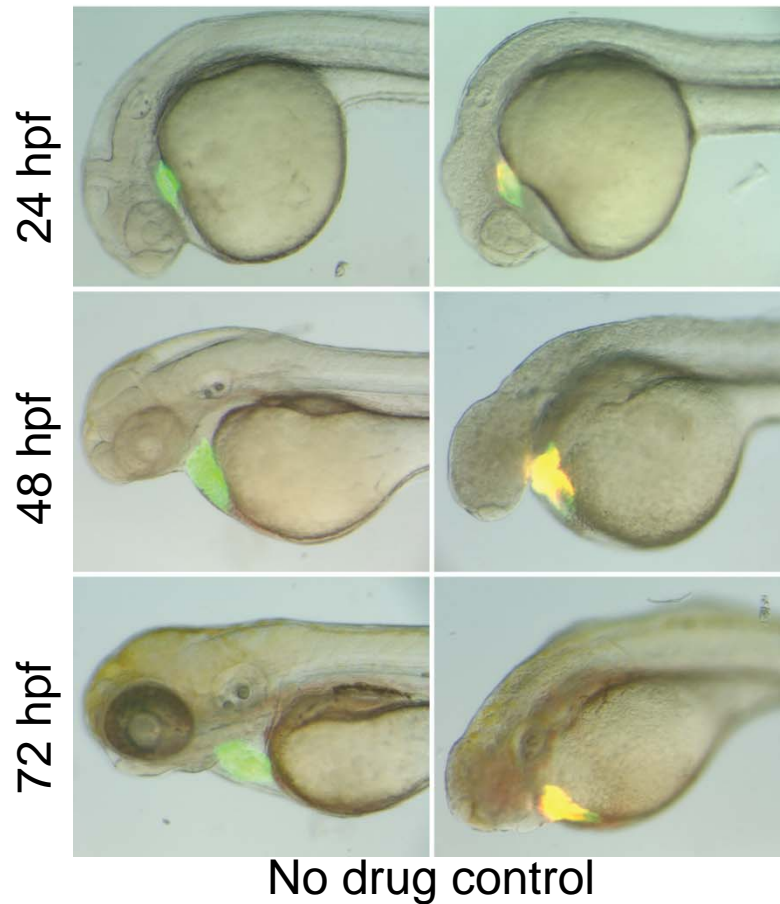
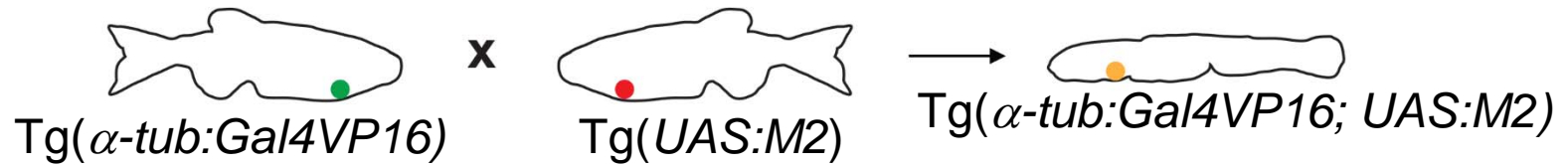
Generating and characterizing Tg(*UAS:M2*) lines



No drug control

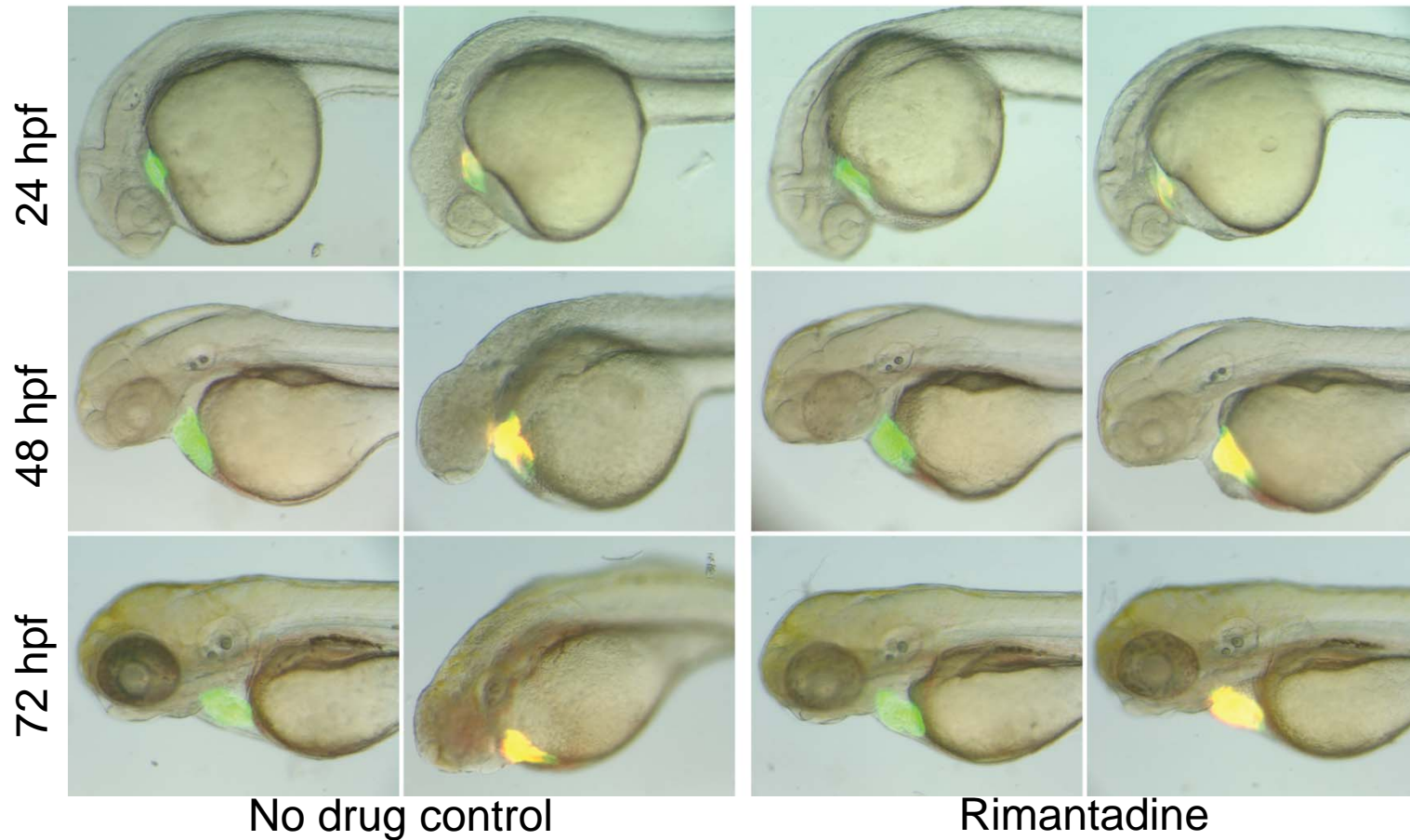
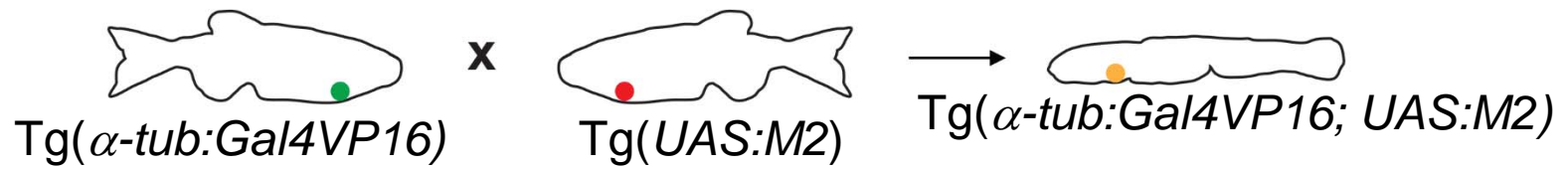
n = 30-50 larvae for each condition

Generating and characterizing Tg(*UAS:M2*) lines



n = 30-50 larvae for each condition

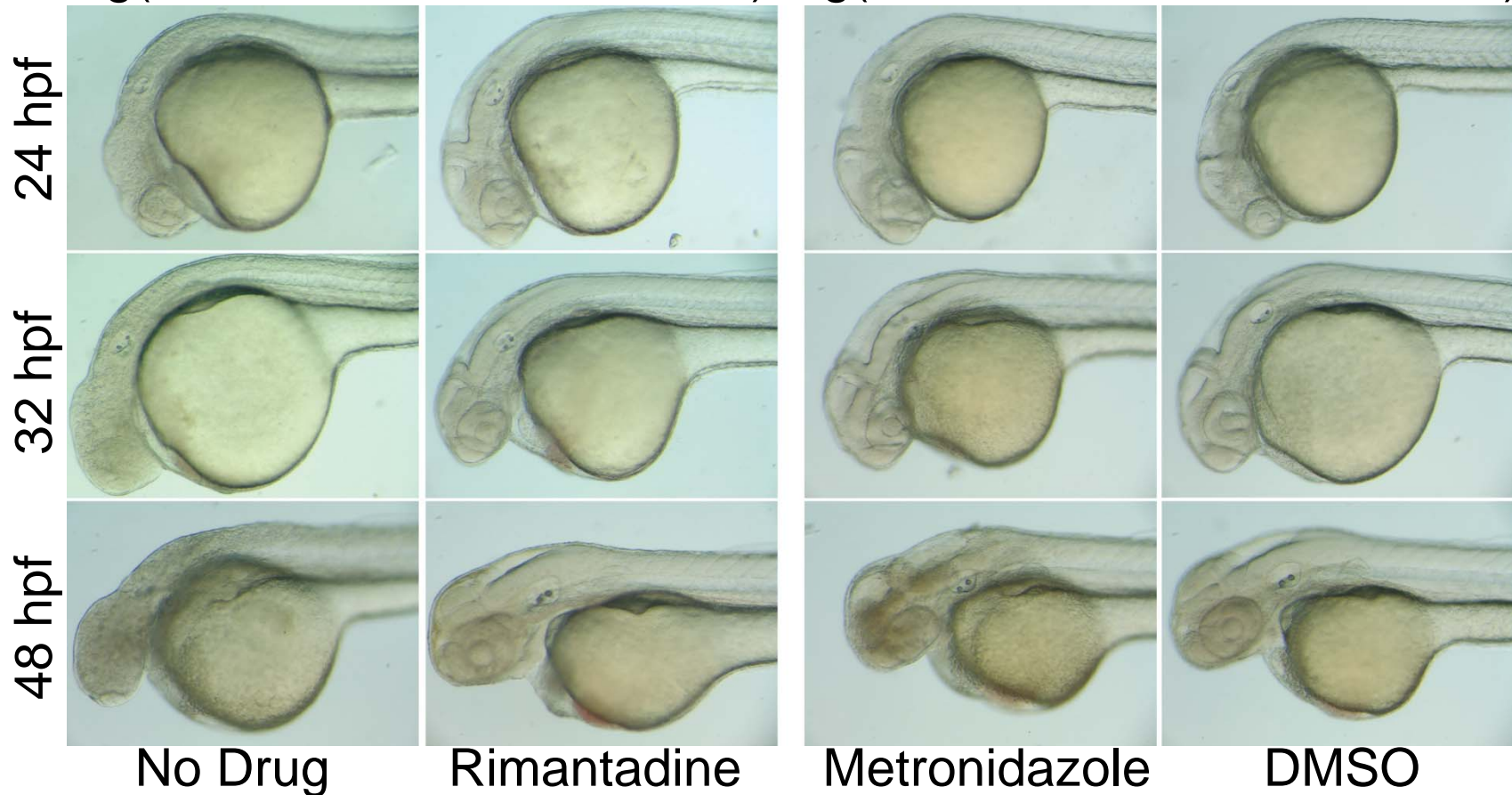
Generating and characterizing Tg(*UAS:M2*) lines



n = 30-50 larvae for each condition

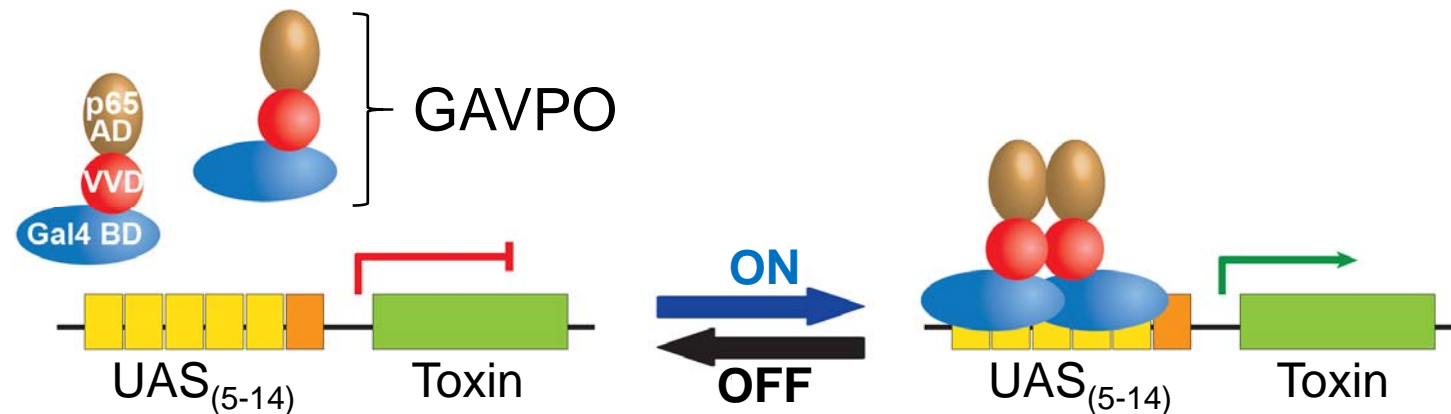
M2-induced defects are more pronounced than NTR-mediated ablation

Tg(α -tub:Gal4VP16; UAS:M2) *Tg(α -tub:Gal4VP16; UAS:NTR)*



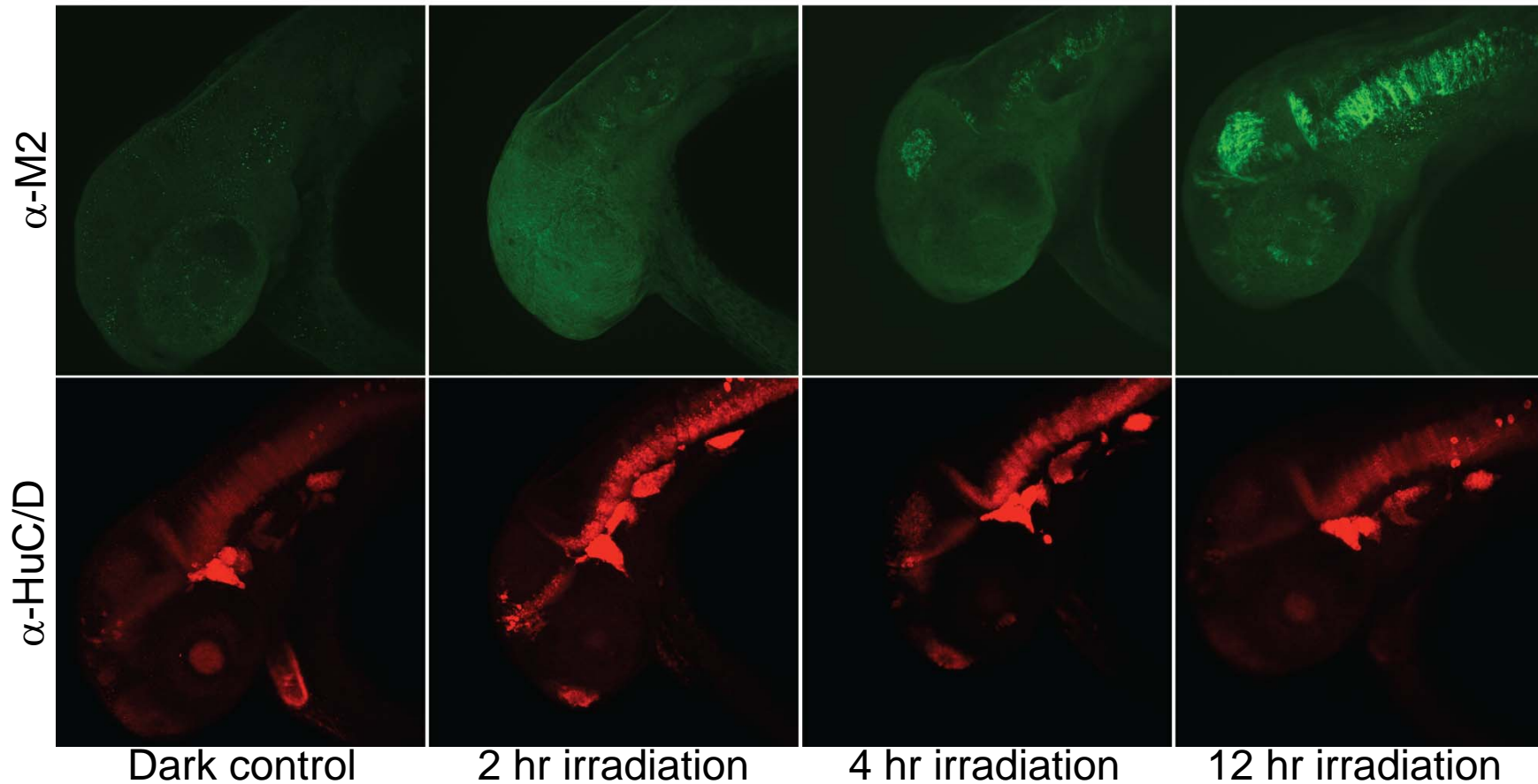
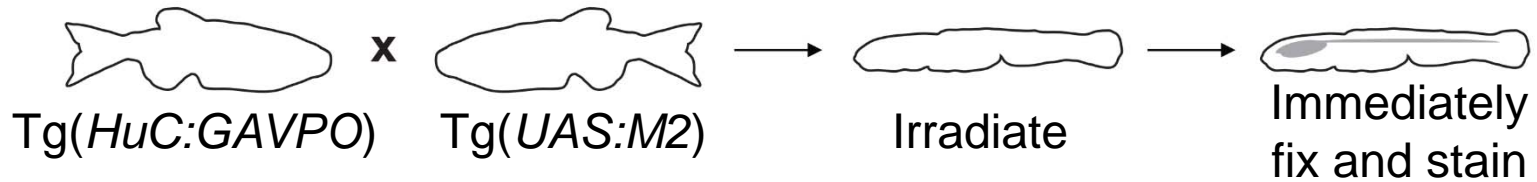
n = 30-50 larvae for each condition

Using light to spatiotemporally ablate CNS cells



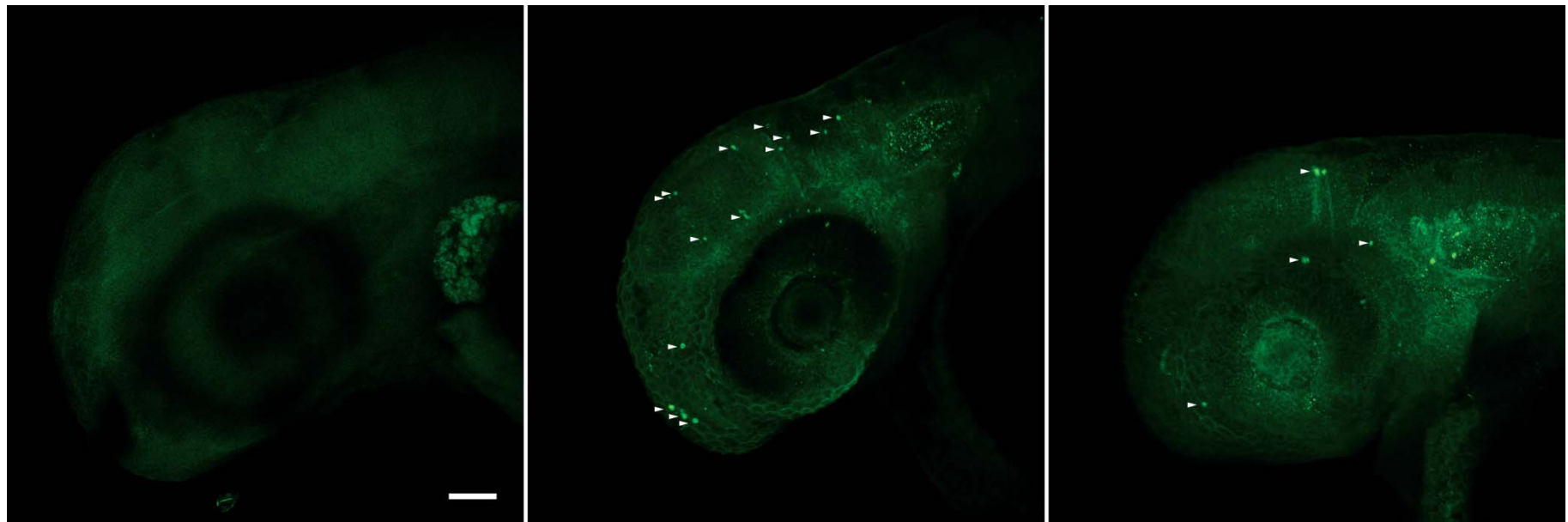
- ✓ Does GAVPO drive transcription in zebrafish?
- ✓ Can we model neural death with toxins?
- ❑ Can we induce neural death using the GAVPO system?

Irradiation induces M2 expression



n = 10-20 larvae for each condition

Irradiation induces activation of Caspase-3 which can be attenuated with rimantadine



Dark control

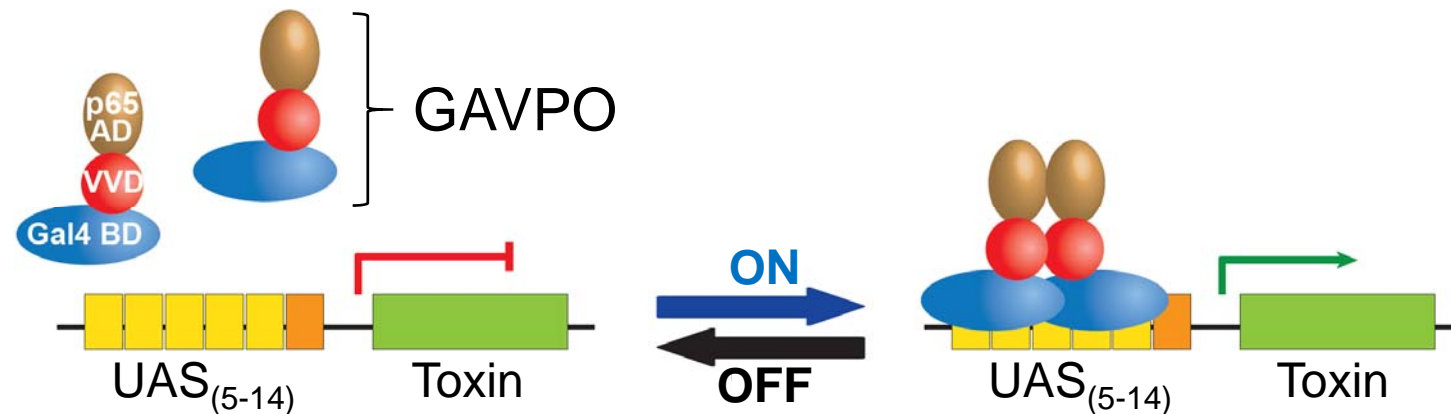
12 hr irradiation

12 hr irradiation
+ rimantadine

Global irradiation (470-nm light 3 mW/cm²)

n = 10-20 larvae for each condition

Using light to spatiotemporally ablate CNS cells



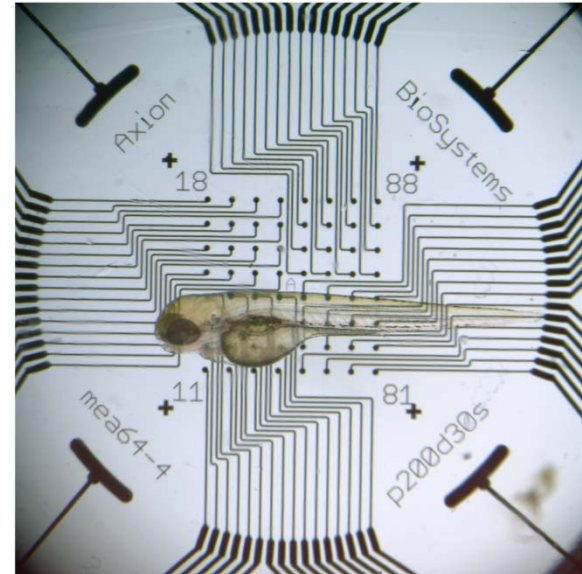
- ✓ Does GAVPO drive transcription in zebrafish?
- ✓ Can we model neural death with toxins?
- ✓ Can we induce neural death using the GAVPO system?

**How does the nervous system response
to optogenetic ablation compare to
manual injury?**

Comparing neural ablation to traditional spinal cord injury in zebrafish



Maestro MEA system
(Axion BioSystems)



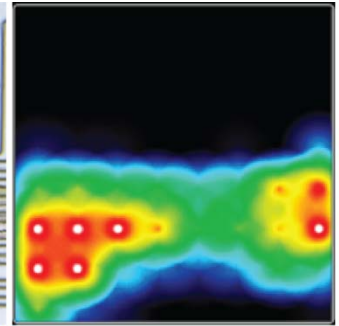
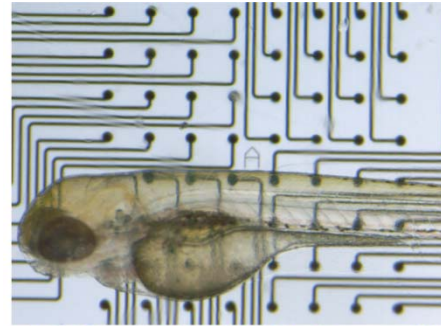
Comparing neural ablation to traditional spinal cord injury in zebrafish



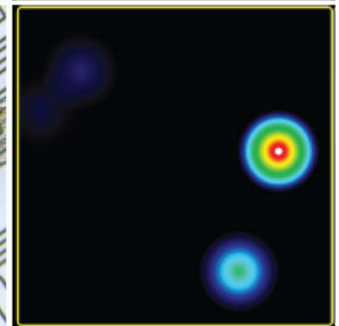
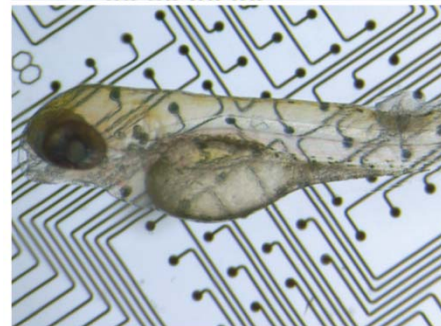
Maestro MEA system
(Axion BioSystems)



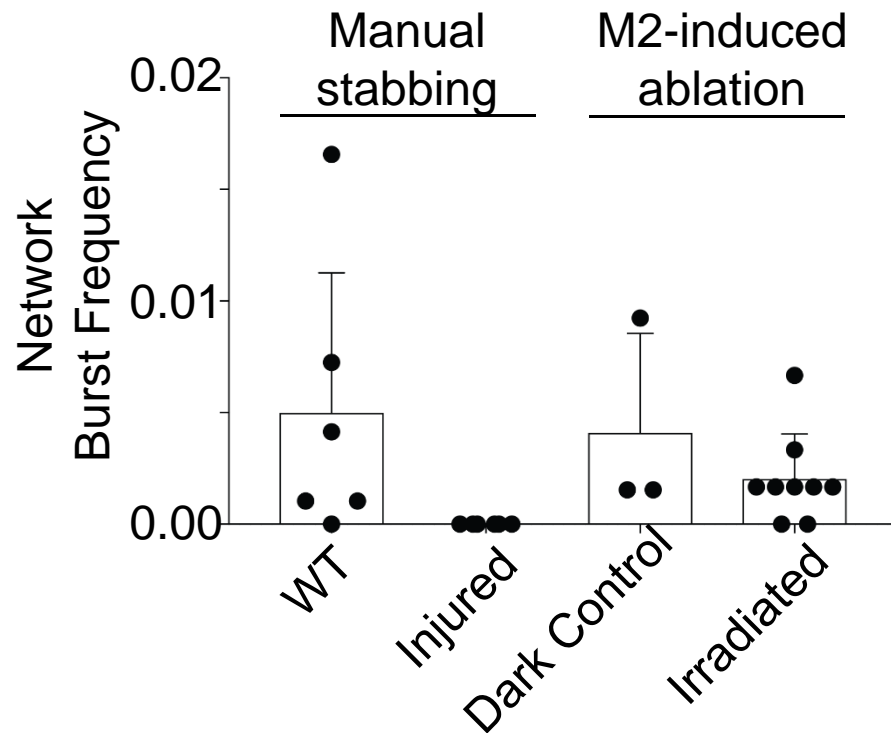
Control



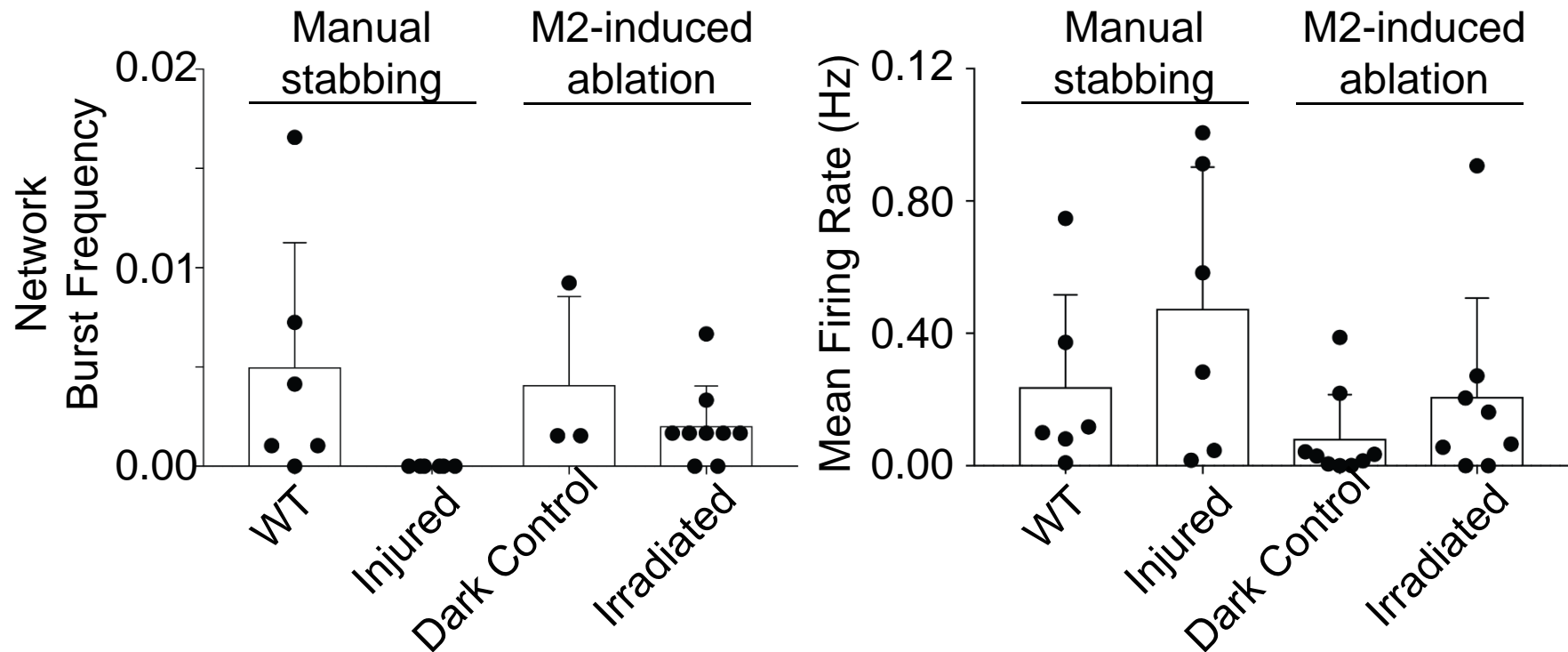
Stabbed



Injury reduces network bursting and increases firing frequency



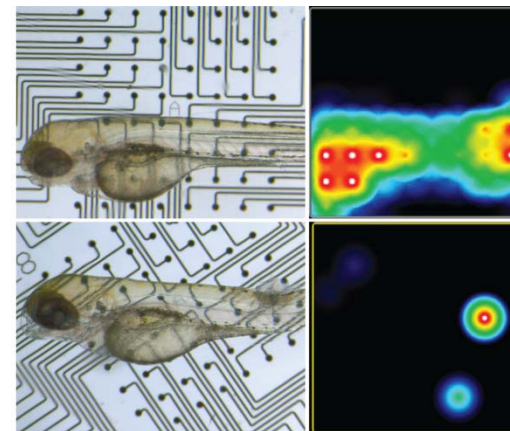
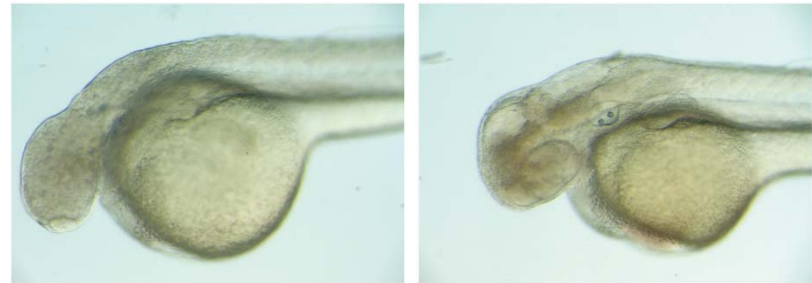
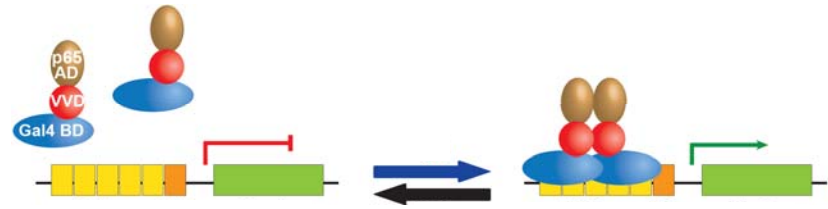
Injury reduces network bursting and increases firing frequency



n = 3-10 larvae for each condition

Summary and future directions

- Developed transgenic zebrafish for light-inducible gene expression
- Used multiple toxins to model secondary injury
- Studying neuronal circuit-connectivity and physiology and SCI pathology



Acknowledgements

- Chen Laboratory
 - **James Chen**
 - Moe Alnaqid
 - Patrick Piza
- Plant Laboratory
 - **Giles Plant**
- Axion Biosystems
 - Stacie Chvatal
 - Emily Matheu

