

Boston Interactive Glial Group (BIGG) Meeting

University of Massachusetts Medical School
Friday, May 17, 2019

Organizer:

– Dorothy Schafer (University of Massachusetts Medical School)

Keynote speakers:

– Dwight Bergles (Johns Hopkins University)

Speaker line-up to be announced.

Topics:

Cell biology of glial cell development

Glial cell-mediated regulation of neuronal structure and physiology

Glial-mediated mechanisms of disease