

## New Fluorescent Tools to Monitor Cell Stress and Toxicity



Time (hours)

Kevin Harlen, Thom Hughes, Anne Marie Quinn Montana Molecular, Bozeman, MT montanamolecular.com

and mutant versions of either the rhodopsin or SOD proteins. The bar plot compares data gathered on a plate reader (n=3 wells). The box plot compares imaging data (n=287 - 359 cells depending on condition).



Fellowship Program NSF Award IIP-1552305.