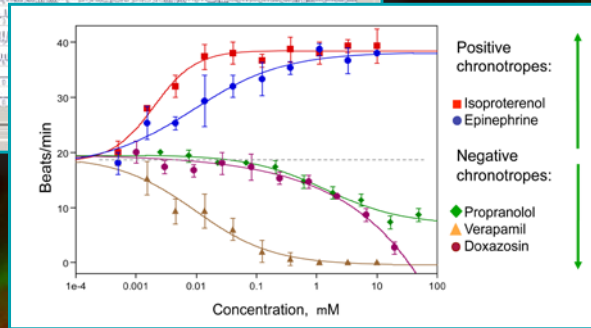


Peak  
Freq = 7.2 bpm  
Rise = 0.53 sec  
Decay = 1.81 sec

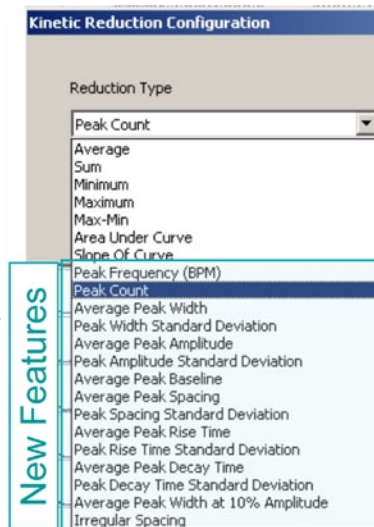


## Introducing ScreenWorks® Peak Pro™ Software for FLIPR® Tetra Cellular Screening Systems

Visualize and analyze the temporal response curves associated with changes in intracellular calcium ion concentration. The new analysis features simplify the characterization of calcium oscillations in primary and stem cells, such as beating cardiomyocytes.

### Enables comprehensive analysis of multi-peak data sets on the FLIPR Tetra System

- Delivers 15 new kinetic analysis parameters to fully characterize data with multiple peaks per measurement
- Measure the impact of pharmacological compounds on the rate and magnitude of beating cardiomyocytes



### Flexibility to optimize peak detection parameters

- Default peak detection parameters easily handle analysis of most cardiomyocyte data
- Configure Peak Detection menu adds ability to further optimize peak detection and capture peak measurements of a variety of size and shape

