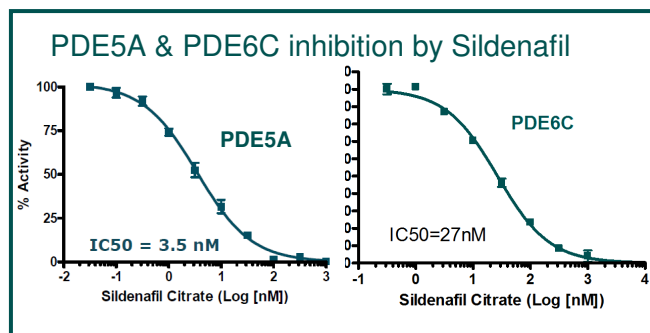
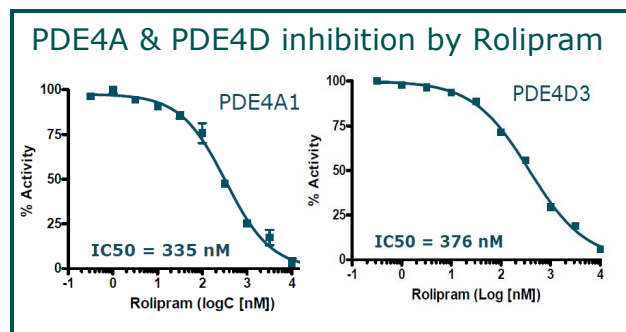
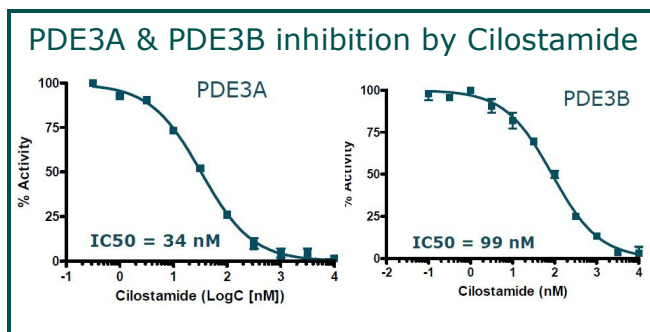
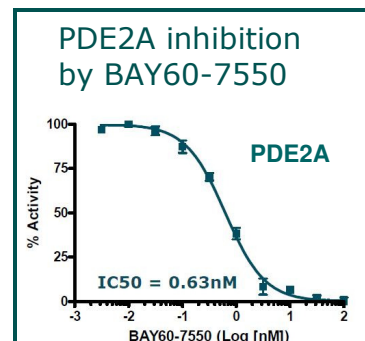
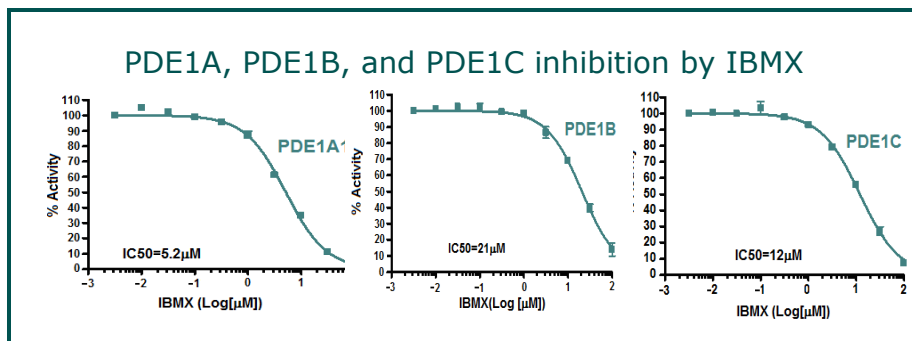


PDE Profiling



Determine IC₅₀ values or screen your compounds against our panel of over 20 different PDE isozymes!

