

Binding kinetics of MK-5108 on AURA

