

Binding kinetics of crizotinib on PYK2

PYK2 Crizotinib

