

## Binding kinetics of crizotinib on PYK2 (PTK2B)

