

Binding kinetics of trametinib on MAP2K1 inactive

