

Binding kinetics of alectinib on ALK

