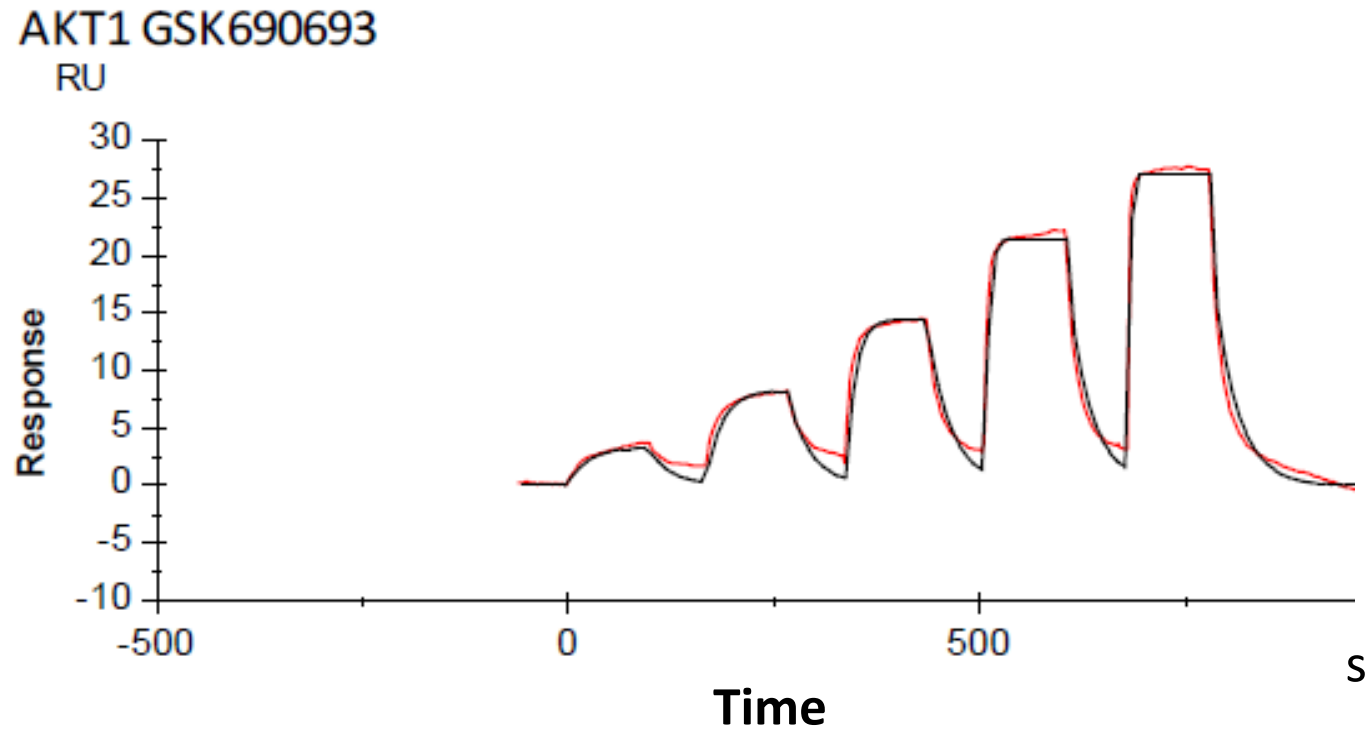


Binding kinetics of GSK690693 on AKT1



The data was generated and provided by
Netherlands Translational Research Center B.V. (NTRC)