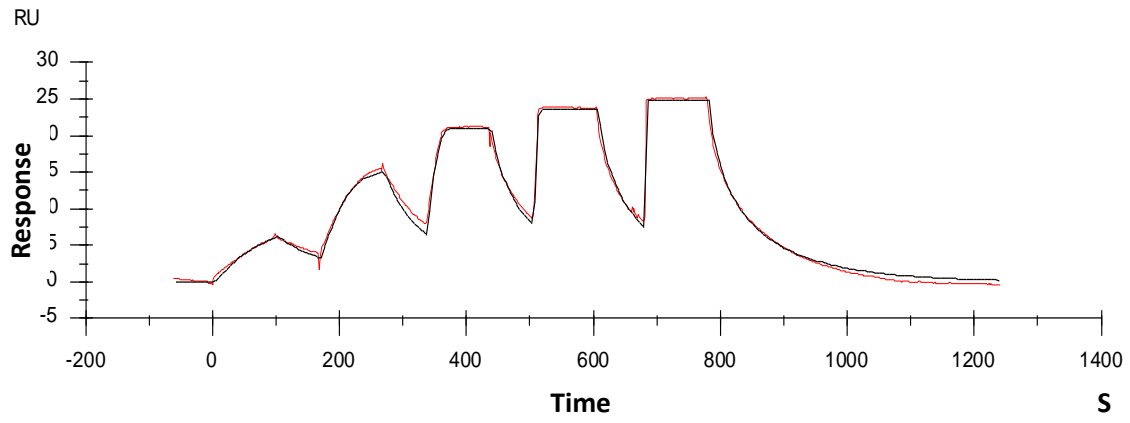


Binding kinetics of A-674563 on BTN-PKAC α (PRKACA)

BTN-PKAC α (PRKACA) A-674563



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