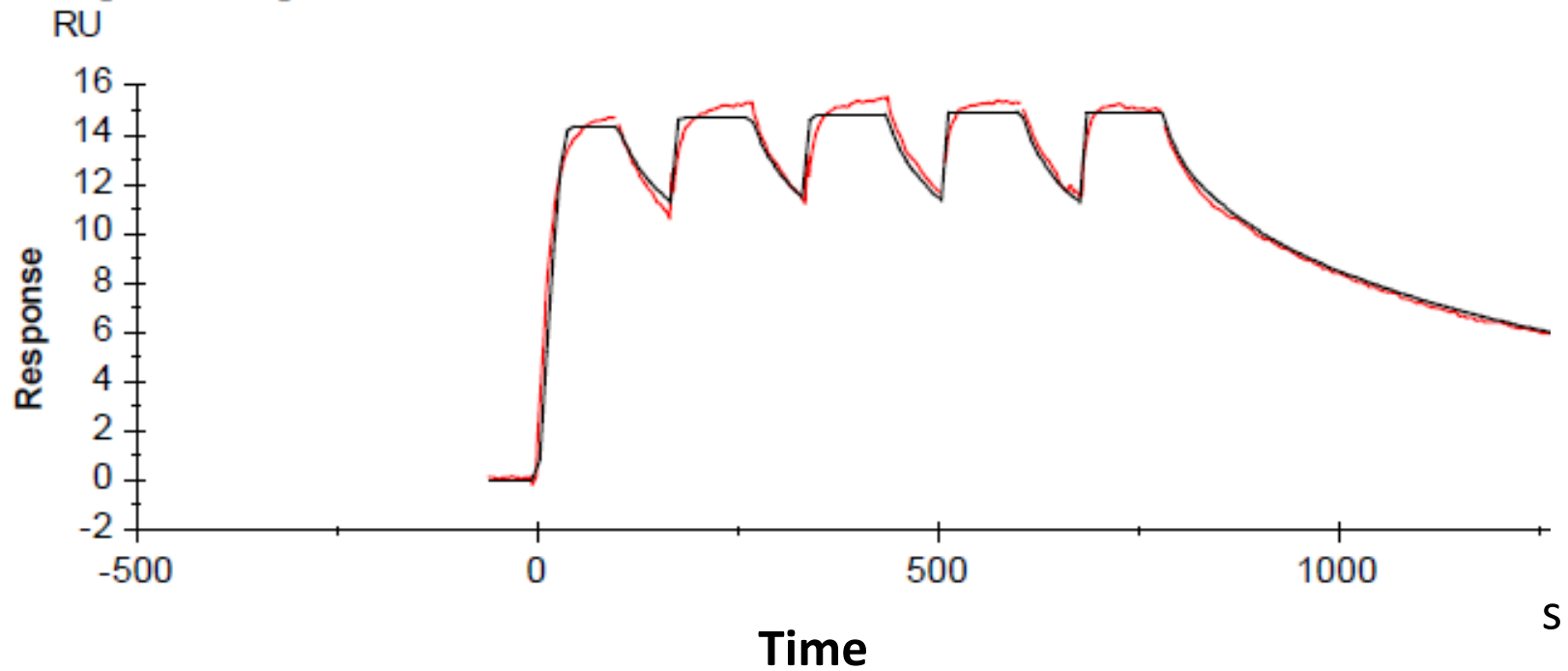


# Binding kinetics of cabozantinib on ALK[L1196M]

## ALK[L1196M] Cabozantinib



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