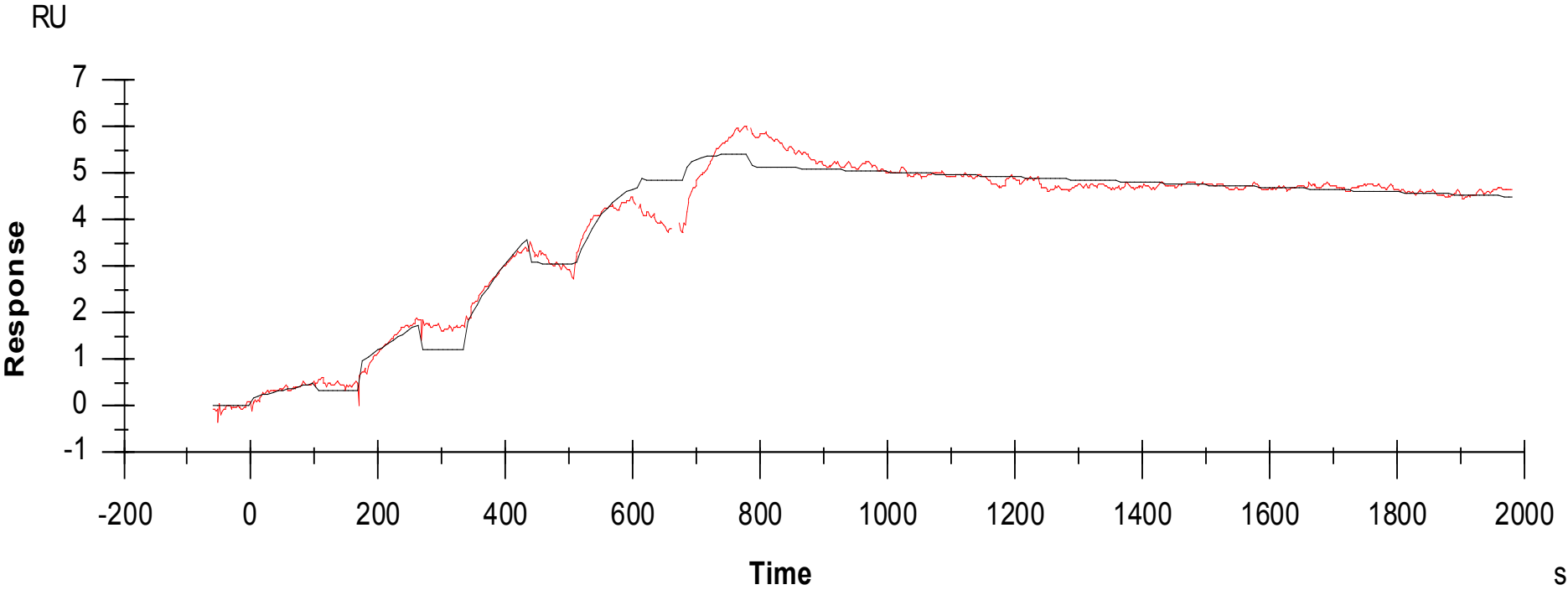


Binding kinetics of AZD1775 on WEE1



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