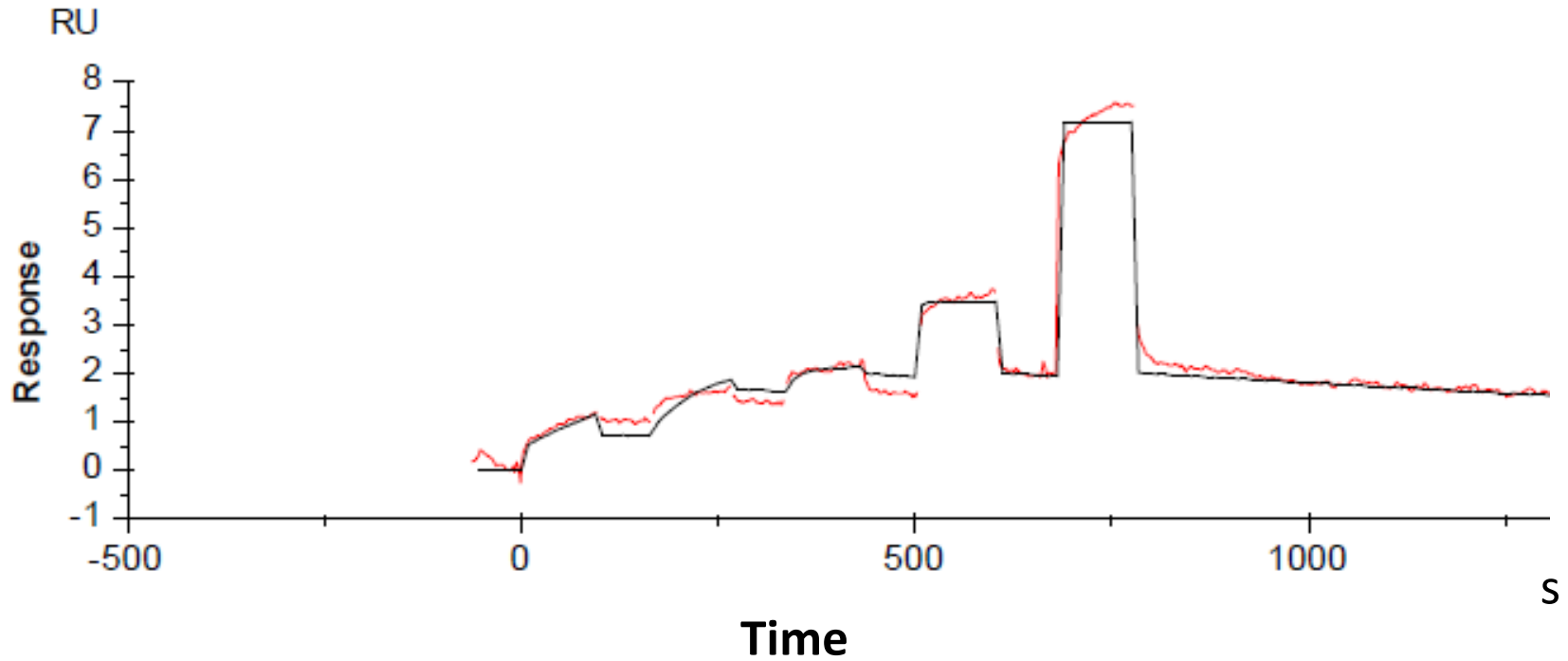


Binding kinetics of BI2536 on PLK1

PLK1 BI2536



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