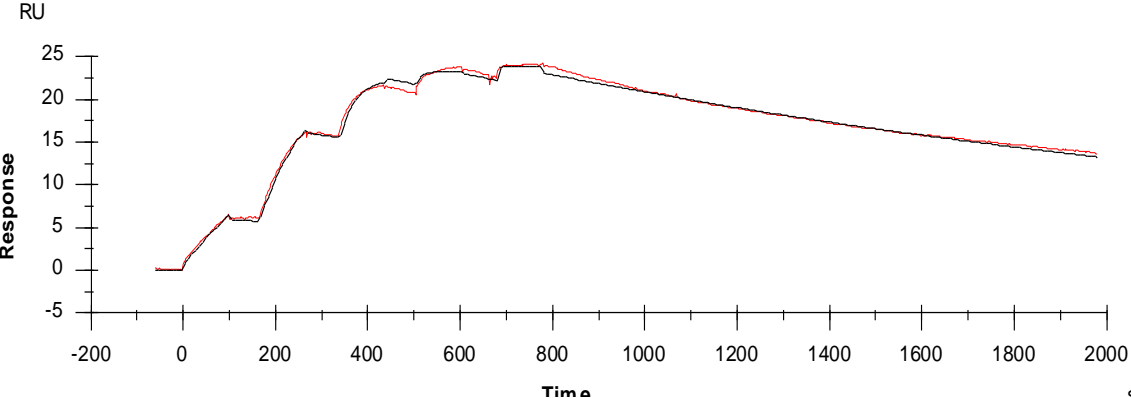


Binding kinetics of BMS-777607 on BTN-MER



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