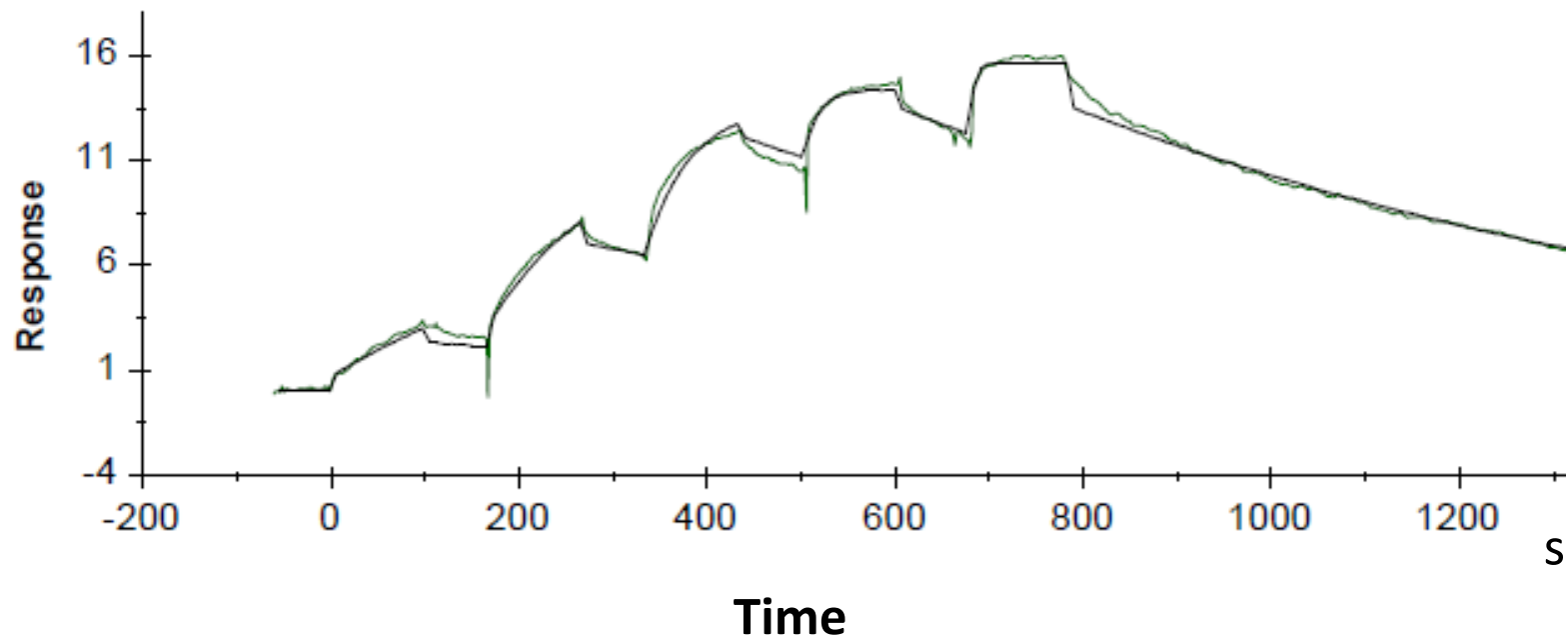


Binding kinetics of vemurafenib on BRAF

vemurafenib on BRAF

RU



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