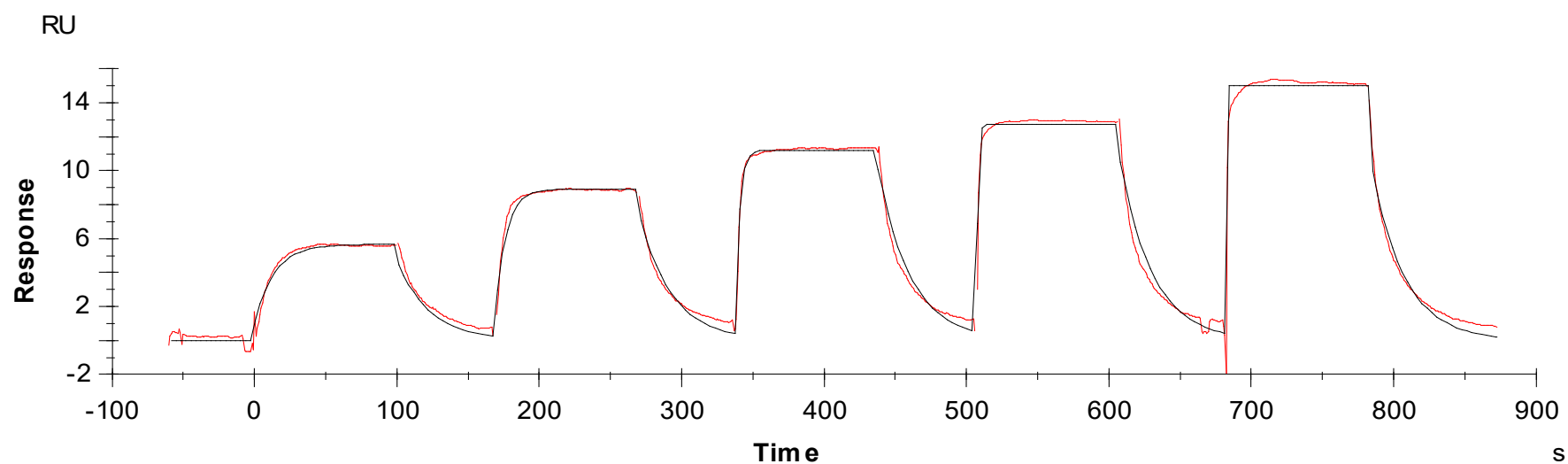
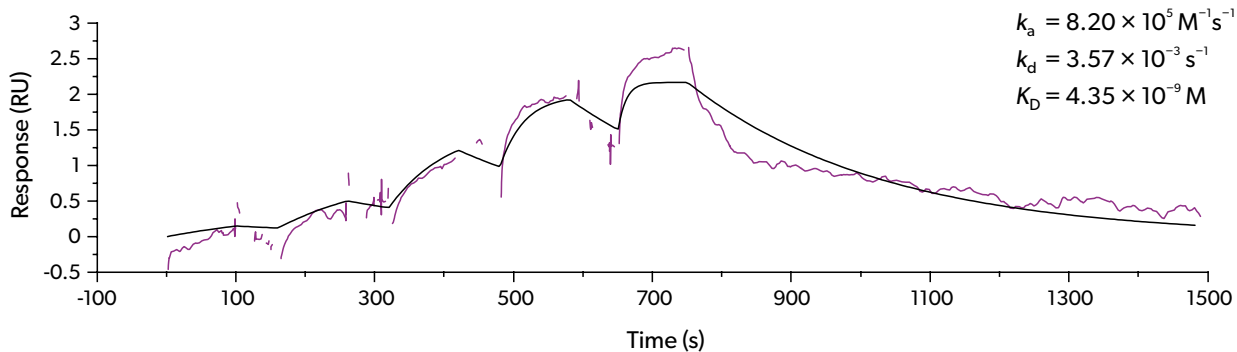


Binding Kinetics of K252A on PKC α (PRKCA)



Data generated and provided by Carna Biosciences, Inc.

Binding kinetics of sotrastaurin on PKC α (PRKCA)



Oncolines

A SYMERES COMPANY

ResidenceTimer™ service

www.oncolines.com

services@oncolines.com