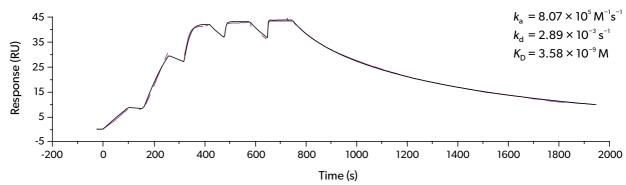
Binding kinetics of sunitinib on FLT3 [non-activated]





ResidenceTimer™ service www.oncolines.com services@oncolines.com