

Binding kinetics of apitolisib on PIK3CB/PIK3R1

