

KinaseSeeker™ - Plasmodium falciparum Panel

Kinase	Organism	GeneDB ID
PfARK1	<i>P. falciparum 3D7</i>	PF3D7_0605300
PfARK3	<i>P. falciparum 3D7</i>	PF3D7_1356800
PfCDPK2	<i>P. falciparum 3D7</i>	PF3D7_0610600
PfCDPK5	<i>P. falciparum 3D7</i>	PF3D7_1337800
PfGSK3	<i>P. falciparum 3D7</i>	PF3D7_0312400
PfNEK1	<i>P. falciparum 3D7</i>	PF3D7_1228300
PfNEK3	<i>P. falciparum 3D7</i>	PF3D7_1201600
PfPK5	<i>P. falciparum 3D7</i>	PF3D7_1356900
PfPK6	<i>P. falciparum 3D7</i>	PF3D7_1337100
PfPK9	<i>P. falciparum 3D7</i>	PF3D7_1337100
PfPKB	<i>P. falciparum 3D7</i>	PF3D7_1246900

To place an order or request a quote,
please contact us at:

Phone: (520) 495-0161
Email: info@luceome.com