

**KinaseSeeker™ - Plasmodium falciparum Panel**

<b>Kinase</b>	<b>Organism</b>	<b>GeneDB ID</b>
PfARK1	<i>P. falciparum 3D7</i>	PF3D7_0605300
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
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](mailto:info@luceome.com)