

Carna's Profiling targets of which reference is to be replaced from January, 2021

Kinase	Current Reference	New Reference
EEF2K	NH125	A-484954
Erk1	5-Iodotubercidin	K252a
Erk1_1mM	5-Iodotubercidin	K252a
Erk2	5-Iodotubercidin	K252a
Erk2_1mM	5-Iodotubercidin	K252a
HPK1	K252b	K252a
JNK1	SP600125	K252a
JNK1_1mM	SP600125	K252a
JNK2	SP600125	K252a
JNK2_1mM	SP600125	K252a
JNK3	SP600125	K252a
JNK3_1mM	SP600125	K252a
MAPKAPK3	K252b	K252a
MINK	K252b	K252a
MINK_1mM	K252b	K252a
MSSK1	Staurosporine	K252a
PLK1	Staurosporine	GW843682X
PLK1_1mM	Staurosporine	GW843682X
PLK2 (IMAP)	K252b	GW843682X
PLK3	K252b	GW843682X
PLK3_1mM	K252b	GW843682X
WNK1	Staurosporine	K252a
WNK2	Staurosporine	K252a
WNK3	Staurosporine	K252a

This change will be reflected on our Profiling Book as of Jan. 1, 2021