

Alignment of Subdomain III Region in Human Cyclin-Dependent Kinases

Subdomain III

CDK1	EEEGVPSTAI REIS LLKELR
CDK2	ETEGVPSTAI REIS LLKELN
CDK3	EMEGVPSTAI REIS LLKELK
CDK4	GGGGLPISTV RE VALLRRLE
CDK5	DDEGVPS SA LI RE ICLLKELK
CDK6	GE EG MP L STI RE VAVLRHLE
CDK7	AKDGINRTAL RE IKLLQELS
CDK8	EGTG IS MSA CRE IAL L RELK