

Phosphorylation Sites in Cyclin-Dependent Kinases

	Subdomain I	Subdomain VIII
Cdc2 (human)	KIGEG ^T YGVV	VY ^T HEVVTLWYRSPEVL
Cdk2 (human)	GIGEG ^T YGVV	TY ^T HEVVTLWYRAPEIL
Cdk3 (human)	KIGEG ^T YGVV	TY ^T HEVVTLWYRAPEIL
Cdk4 (mouse)	EIGVGAYGTV	AL ^T PVVTLWYRAPEVL
Cdk5 (human)	KIGEG ^T YGTV	CY ^S AEVVTLWYRPPDVL
Cdk6 (human)	KIGEG ^S YGVV	YY ^T TDYVATRWYRSPELL
Cdk7 (mouse)	FLGRGQFATV	AY ^T HQVVTRWYRAPELL

T14 Inhibitory phosphorylation

Y15 **K71**



Stimulatory

T161 phosphorylation



ATP-binding
domain

Protein kinase
activation loop