

Anti-Cyclin G1

(Catalog # A-0010)

Background:	Cyclin G contains a typical cyclin box at its N-terminus but no destruction box. Cyclin G levels rise when DNA is damaged. Cyclin G seems to promote apoptosis, although its function is not exactly known. Cyclin G is noted to regulate the levels of p53 and to be involved in tumorigenesis.
Concentration:	0.2 mg/ml
Description:	Rabbit polyclonal antibody to cyclin G1. Epitope corresponding to 46 amino acids of the N terminal region of cyclin G1.
Purification:	Protein A Purified
Specificity:	Mouse and human
Isotype:	IgG
Formulation:	0.05% Sodium Azide, 0.1% Gelatin, PBS
Storage:	4 °C, stable for 1 year from the date of shipment.
Application:	Western blot: 1:200-1:1000 Immunofluorescence: 1:100-1: 500 Immunohistochemistry: 1:100-1:500 ELISA: 1:1000-1:2000 Immunoprecipitation: 2 µg/10 ⁶ cells
Research Use:	Research use only.

Ordering Information

Products	Size	Cat. No.
Anti-Cyclin G1	50 µg	A-0010-050
	100 µg	A-0010-100

Available Related Products	Size	Cat. No.
Anti-Cyclin G2	50 µg	A-0011-050
	100 µg	A-0011-050

This product is for research purposes only. Not intended for use in diagnostic procedures.