

Anti-Cyclin G2

(Catalog # A-0011)

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:	0.5 mg/ml
Description:	Goat polyclonal antibody to cyclin G2.
Purification:	Protein A Purified
Specificity:	Mouse and human
Isotype:	IgG
MW:	39 KDa
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 ELISA: 1:1000-1: 2000
Research Use:	Research use only.

Ordering Information

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

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

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