







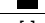
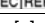
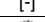
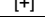



**Rabbit (polyclonal)
Anti-Murine p21^{WAF1}/Cip1/Sdi1/Pic1/CAP20
(Mitotic Inhibitor)
Unconjugated
PRODUCT ANALYSIS SHEET**

Catalog Number:	AMZ0402
Lot Number:	See product label
Expiration Date:	See product label.
Quantity/Volume:	500 µg/0.5 mL
Form of Antibody:	Purified immunoglobulin in phosphate buffered saline, pH 7.4, with 0.2% bovine serum albumin.
Preservation:	0.09% sodium azide (Caution: sodium azide is a poisonous and hazardous substance. Handle with care and dispose of properly.)
Purification:	Purified from rabbit serum by Protein A chromatography.
Immunogen:	A synthetic peptide corresponding to aa 2-14 (Cys-SNPGDVRPVPHRSK) from the N-terminus of mouse p21 ^{WAF1} . The peptide was coupled to KLH.
Specificity:	This antibody recognizes a protein of 21 kDa, identified as the p21 ^{WAF1} /Cip1/Sdi1/Pic1 tumor suppressor protein, a tumor suppressor involved in the pathogenesis of a variety of malignancies. AMZ0402 is highly specific to p21 and shows no cross-reactivity to other closely related mitotic inhibitors. Expression of p21 ^{WAF1} is induced by wild type, but not mutant, p53 suppressor protein. The p21 ^{WAF1} protein binds to cyclin/cdk complexes and inhibits their kinase activity, arresting cell cycle progression. It also binds to PCNA (proliferating cell nuclear antigen) and blocks DNA replication but not the DNA repair process. Inactivation of p53 is believed to disrupt the p53-dependent expression of p21 ^{WAF1} .
Species Reactivity:	Mouse and rat p21 ^{WAF1} ; does not cross-react with human. Other species were not tested.
Applications:	This antibody is suitable for use in kinase assays, immunoprecipitation (co-precipitation of cdk4/cdk6) and immunohistochemistry with cryostat and formalin-fixed paraffin embedded tissues. Use of this antibody with formalin-fixed paraffin embedded tissues requires boiling the tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 minutes, followed by cooling for 20 minutes at room temperature.
Suggested Working Dilutions:	For immunoprecipitation, use 10 µL/mg of protein lysate. The optimal antibody concentration should be determined for each specific application.
Recommended Positive Control:	3T3-L1 cells; rat or mouse embryonic fibroblast cells treated with DNA damaging agents; skin, colon or breast carcinoma.
Storage:	Store at 2-8°C.

Explanation of symbols

Symbol	Description	Symbol	Description
	Catalogue Number		Batch code
	Research Use Only		In vitro diagnostic medical device
	Use by		Temperature limitation
	Manufacturer		European Community authorised representative
	Without, does not contain		With, contains
	Protect from light		Consult accompanying documents
	Directs the user to consult instructions for use (IFU), accompanying the product.		

For Research Use Only. CAUTION: Not for human or animal therapeutic or diagnostic use.

www.invitrogen.com

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288 • E-mail: techsupport@invitrogen.com

PI AMZ0402

(Rev 04/10) DCC-10-0604

Important Licensing Information - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, www.invitrogen.com). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.