

RAD23A Antibody [#C13113]

Catalog Number:	C13113
Concentration:	1mg/ml
Swiss-Prot No.:	P54725
Other Names:	UV excision repair protein RAD23 homolog A, hHR23A, RAD23A
Storage/Stability:	Store at -20°C/1 year
Form of Antibody:	Rabbit IgG in phosphate buffered saline (without Mg ²⁺ and Ca ²⁺), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
Immunogen:	The antiserum was produced against synthesized peptide derived from internal of human RAD23A.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	RAD23A antibody detects endogenous levels of total RAD23A protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Masutani C., EMBO J. 13:1831-1843(1994). Grimwood J., Nature 428:529-535(2004). Withers-Ward E.S., J. Virol. 71:9732-9742(1997).