

## Bloom Syndrome Antibody [#B0778]

<b>Catalog Number:</b>	B0778
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P54132
<b>Other Names:</b>	Bloom's syndrome protein, DNA helicase, RecQ-like, type 2, EC 3.6.1.-, RECQ2, RECQL3, RecQ protein-like 3
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from Internal of human Bloom Syndrome.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	Bloom Syndrome antibody detects endogenous levels of total Bloom Syndrome protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%)
<b>Applications:</b>	IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:5000
<b>References:</b>	V. Ashutosh Rao. AACR Meeting Abstracts, Mar 2004; 2004: 1244. Xiaobing Du. Mol. Cell. Biol., Oct 2004; 24: 8437 - 8446. Patricia L. Opresko. J. Biol. Chem., Oct 2002; 277: 41110 - 41119.