

MITF (Ab-180/73) Antibody [#B0512]

Catalog Number:	B0512
Concentration:	1mg/ml
Swiss-Prot No.:	O75030
Other Names:	MITF M1, Microphthalmia-associated transcription factor
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 non-phosphopeptide derived from human MITF around the phosphorylation site of serine 180/73 (P-N-S ^P -P-M).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MITF (Ab-180/73) antibody detects endogenous levels of total MITF protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:10000
References:	Jón H. Hallsson, Genetics, May 2004; 167: 233 - 241. Eiichi Morii, Blood, Apr 2001; 97: 2038 - 2044. Eiichi Morii, Blood, Sep 2004; 104: 1656 - 1661.