

IF3EI Antibody [#C11925]

Catalog Number:	C11925
Concentration:	1mg/ml
Swiss-Prot No.:	Q9Y262
Other Names:	EIF3S6IP, Eukaryotic translation initiation factor 3 subunit 6 interacting protein, HSPC021/HSPC025, IF3I
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 N-terminal of human IF3EI.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	IF3EI antibody detects endogenous levels of total IF3EI protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Zhang Q.-H., Genome Res. 10:1546-1560(2000). Collins J.E., Genome Biol. 5:RESEARCH84.1-RESEARCH84.11(2004). Morris-Desbois C., J. Biol. Chem. 276:45988-45995(2001).