

LRRC41 Antibody [#C16455]

Catalog Number:	C16455
Concentration:	1mg/ml
Swiss-Prot No.:	Q15345
Other Names:	Leucine-rich repeat-containing protein 41, Protein Muf1
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 LRRC41.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	LRRC41 antibody detects endogenous levels of total LRRC41 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Gregory S.G., Nature 441:315-321(2006). The MGC Project Team, Genome Res. 14:2121-2127(2004). Bechtel S., BMC Genomics 8:399-399(2007).