

TRAPPC3 Antibody [#C19146]

Catalog Number:	C19146
Concentration:	1mg/ml
Swiss-Prot No.:	O43617
Other Names:	Trafficking protein particle complex subunit 3; BET3 homolog; TRAPPC3; BET3; CDABP0066
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 TRAPPC3.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	TRAPPC3 antibody detects endogenous levels of total TRAPPC3 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Sacher M., EMBO J. 17:2494-2503(1998). Eva L., Submitted (JAN-1998) to the EMBL/GenBank/DDBJ databases. Zhou J., Submitted (JUL-2000) to the EMBL/GenBank/DDBJ databases.