

## DNAL1 Antibody [#C15512]

**Catalog Number:** C15512

**Concentration:** 1mg/ml

**Swiss-Prot No.:** Q4LDG9

**Other Names:** C14orf168; dynein light chain 1, axonemal

**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 DNAL1.

**Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.

**Specificity:** DNAL1 antibody detects endogenous levels of total DNAL1 protein.

**Reactivity:** Human, Mouse, Rat

**Applications:** WB: 1:500~1:1000 ELISA: 1:40000

**References:** Horvath J., Am. J. Respir. Cell Mol. Biol. 33:41-47(2005).

The German cDNA consortium; Submitted (JAN-2005) to the EMBL/GenBank/DDBJ databases.

The MGC Project Team; Genome Res. 14:2121-2127(2004).