

Dysferlin Antibody [#C0176]

- Catalog Number:** C0176
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- Swiss-Prot No. :** O75923
- All Names:** dysferlin, limb girdle muscular dystrophy 2B (autosomal recessive)
- Form of Antibody:** Rabbit IgG in phosphate buffered saline (without Mg²⁺ and Ca²⁺), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
- Storage/Stability:** Store at -20°C/1 year
- Immunogen:** The antiserum was produced against synthesized peptide derived from human Dysferlin.
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity** Dysferlin antibody detects endogenous levels of total Dysferlin protein.
- Reactivity:** Human, Mouse
- Applications:** WB: 1:300~1:1000 ELISA: 1:10000
- References:** Delia J. Hernández-Deviez, Hum. Mol. Genet., Jan 2006; 15: 129 - 142.
Yanchao Huang, FASEB J, Mar 2007; 21: 732 - 742.
Mengfatt Ho, Hum. Mol. Genet., Sep 2004; 13: 1999 - 2010.
Eriko Fujita, Hum. Mol. Genet., Mar 2007; 16: 618 - 629.