

## ABCB7 Antibody [#C14621]

|                           |   |
|---------------------------|---|
| <b>Catalog Number:</b>    | C14621  |
| <b>Concentration:</b>     | 1mg/ml  |
| <b>Swiss-Prot No.:</b>    | O75027  |
| <b>Other Names:</b>       | ABC transporter 7 protein; ABC7; Anemia, sideroblastic, with spinocerebellar ataxia; ASAT; Atm1p; ATP-binding cassette 7; ATP-binding cassette sub-family B member 7, mitochondrial; ATP-binding cassette transporter 7; ATP-binding cassette, sub-family B (MDR/TAP), member 7 |
| <b>Storage/Stability:</b> | Store at -20°C/1 year   |
| <b>Form of Antibody:</b>  | Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.  |
| <b>Immunogen:</b>         | The antiserum was produced against synthesized peptide derived from internal of human ABCB7.  |
| <b>Purification:</b>      | The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.   |
| <b>Specificity:</b>       | ABCB7 antibody detects endogenous levels of total ABCB7 protein.  |
| <b>Reactivity:</b>        | Human (Identities = 100%, Positives = 100%)   |
| <b>Applications:</b>      | IHC: 1:50~1:100 ELISA: 1:40000  |
| <b>References:</b>        | Bekri S., Blood 96:3256-3264(2000).<br>Ross M.T., Nature 434:325-337(2005).<br>Maguire A., Br. J. Haematol. 115:910-917(2001).  |