

## FANCD2 (Ab-222) [#B0427]

- Catalog Number:** B0427-1, B0427-2, B0427
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- Swiss-Prot No. :** Q9BXW9
- All Names:** FACD2, FANCD2, Fanconi anemia complementation group D2 protein
- Form of Antibody:** 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.
- Storage/Stability:** Store at -20°C/1 year
- Immunogen:** The antiserum was produced against synthesized non-phosphopeptide derived from human FANCD2 around the phosphorylation site of serine 222 (G-D-S-Q-H).
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity** FANCD2 (Ab-222) antibody detects endogenous levels of total FANCD2 protein.
- Reactivity:** Human
- Applications:** WB: 1:500~1:1000    IHC: 1:50~1:100  
ELISA: 1:10000
- References:** Shobbir Hussain, Hum. Mol. Genet., Jun 2004; 13: 1241 - 1248.  
Rocio Montes de Oca, Blood, Feb 2005; 105: 1003 - 1009.  
XiaoZhe Wang, Mol. Cell. Biol., Jul 2004; 24: 5850 - 5862.  
Gary P. H. Ho, Mol. Cell. Biol., Sep 2006; 26: 7005 - 7015.