

NRF2 Antibody [#C0279]

Catalog Number: C0279-1, C0279-2, C0279

Amount: 50µg/50µl, 100µg/100µl, 200µg/200µl

Swiss-Prot No. : Q16236

All Names: HEBP1, NF-E2 related factor 2, NF2L2, NFE2-related factor 2, NFE2L2, NFL2_HUMAN, NRF-2, Nuclear factor erythroid 2 related factor 2, Nuclear factor erythroid derived 2 like 2

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 NRF2.

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

Specificity/Sensitivity NRF2 antibody detects endogenous levels of total NRF2 protein.

Reactivity: Human, Mouse, Rat

Applications: WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:5000

References: Rajesh K. Thimmulappa, J. Clin. Invest., Apr 2006; 116: 984 - 995.

Sara B. Cullinan, Mol. Cell. Biol., Oct 2003; 23: 7198 - 7209.

Tirumalai Rangasamy, J. Exp. Med., Jul 2005; 202: 47 - 59.

Wenge Li, J. Biol. Chem., Aug 2005; 280: 28430 - 28438.