

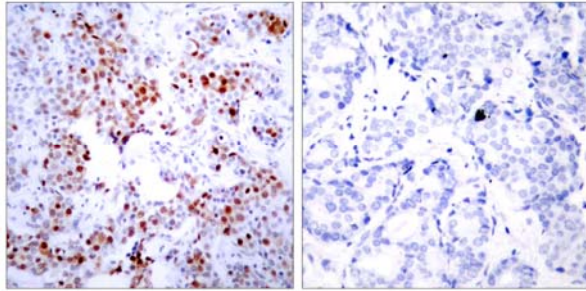


ATF-2 (Ab-62 or 44) Antibody

Catalog Number 21029

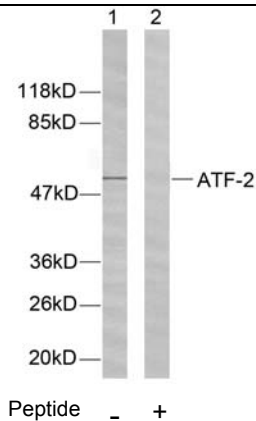
(For Research Use Only)

- Catalog Number:** 21029-1, 21029-2, 21029
Amount: 50µg/50µl, 100µg/100µl, 200µg/200µl
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 non-phosphopeptide derived from human ATF-2 around the phosphorylation site of serine 62 or 44.
Purification: The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity/Sensitivity: ATF-2 (Ab-62 or 44) antibody detects endogenous levels of total ATF-2 protein.
Reactivity: Human, Mouse, Rat
Applications: WB: 1:500~1:1000 IHC: 1:50~1:100
ELISA: 1:10000 IP: Various Dilution
Swiss-Prot No. : P15336
References: Sevilla A, et al. (2004) J Biol Chem. 279(26):27458-27465.
Sakurai A, et al. (1991) Biochem Biophys Res Commun. 181(2): 629-635.
Abdel-Hafiz H A, et al. (1992) Mol Endocrinol. 6: 2079-2089.
Gupta S, et al. (1995) Science. 267: 389-393.
Van Dam H, et al. (1995) EMBO J. 14(8): 1798-1811.



Peptide - +

Immunohistochemical analysis of paraffin-embedded human breast carcinoma tissue using ATF-2 (Ab-62 or 44) antibody (#21029).



Western blot analysis of extracts from HeLa cells using ATF-2 (Ab-62 or 44) antibody (#21029).