

## ARNT Antibody [#C10491]

|                           |   |
|---------------------------|---|
| <b>Catalog Number:</b>    | C10491  |
| <b>Concentration:</b>     | 1mg/ml  |
| <b>Swiss-Prot No.:</b>    | P27540  |
| <b>Other Names:</b>       | Aryl hydrocarbon receptor nuclear translocator, Dioxin receptor, nuclear translocator, HIF-1 beta, Hypoxia-inducible factor 1 beta                                  |
| <b>Storage/Stability:</b> | Store at -20°C/1 year   |
| <b>Form of Antibody:</b>  | Rabbit IgG in phosphate buffered saline (without $Mg^{2+}$ and $Ca^{2+}$ ), 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 ARNT.   |
| <b>Purification:</b>      | The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.   |
| <b>Specificity:</b>       | ARNT antibody detects endogenous levels of total ARNT protein.  |
| <b>Reactivity:</b>        | Human, Mouse, Rat   |
| <b>Applications:</b>      | WB: 1:500~1:1000   ELISA: 1:20000   |
| <b>References:</b>        | Hoffman E.C., Science 252:954-958(1991).<br>Wang G.L., Proc. Natl. Acad. Sci. U.S.A. 92:5510-5514(1995).<br>Nguyen T.A., Arch. Biochem. Biophys. 367:250-257(1999). |