

PGC1- β antibody (pAb)**Catalog Nos:** 39859, 39860**RRID:** AB_2793372**Isotype:** IgG**Application(s):** WB**Reactivity:** Human**Volumes:** 100 μ l, 10 μ l**Purification:** Affinity Purified**Host:** Rabbit**Molecular Weight:** 150 kDa

Background: PGC1- β (Peroxisome proliferator-activated receptor gamma coactivator 1-beta, PPARGC1B, PERC) is a co-activator of transcription by nuclear receptors and transcription factors. PGC1- β participates in modulating the transcriptional activity of PPAR alpha and gamma, ER alpha, NRF1 and GR. It is likely involved in fat oxidation, non-oxidative glucose metabolism and in the regulation of energy expenditure. Expression of PGC1- β is repressed by saturated fatty acids in skeletal muscle cells and induced by insulin. PGC1- β is down-regulated in type 2 diabetes as well as in pre-diabetics. Allelic variation of PGC1- β may contribute to the pathogenesis of obesity, with the widespread Ala-203 variation being a significant risk factor for the development of this diabetes.

Immunogen: This PGC1- β antibody was raised against a peptide within the N-terminal region of human PGC1- β .

Buffer: Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 mM glycine, 0.07 mM EDTA, 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

Applications Validated by Active Motif:

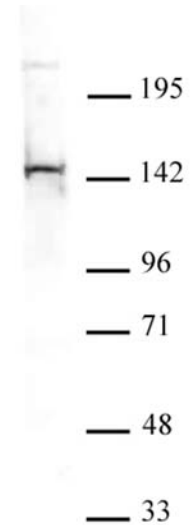
WB*: 1:500 - 1:2,000 dilution

The addition of 0.05% Tween 20 in the blocking buffer and primary antibody incubation buffer is recommended to aid in detection by Western blot. Individual optimization may be required.

*Note: many chromatin-bound proteins are not soluble in a low salt nuclear extract and fractionate to the pellet. Therefore, we recommend a High Salt / Sonication Protocol when preparing nuclear extracts for Western blot.

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.

**PGC1- β antibody tested by Western blot.**

Western blot of K-562 nuclear extract (40 μ g/lane) probed with PGC1- β antibody (pAb) (1:500 dilution).