

Product Information

Description:	Recombinant human PPARD (peroxisome proliferator-activated receptor delta), variant 1, encompassing amino acids 2-441(end). This construct contains an N-terminal His-tag (6xHis). This protein was affinity purified.
Background:	PPARD (peroxisome proliferator-activated receptor delta), also known as NUC1 (nuclear hormone receptor 1), is a hormone receptor of the PPAR group of nuclear receptor superfamily of proteins. It is a ligand-activated transcription factor, present in the nucleus of many cell types, such as heart, lung, brain, and liver. It participates in cellular metabolism homeostasis, being involved in fatty acid uptake, transport and oxidation, insulin secretion and response. Dysfunction of these functions can result in obesity, dyslipidemia, type 2 diabetes, hepatosteatosi and atherosclerosis. All these disorders represent a large burden on both the human body and the health system of industrialized countries, and large efforts have been put into developing clinical strategies to ameliorate them. The development of PPARD agonist was initially seen as a promising therapeutic strategy, however it was found that cancer cells, which use large amounts of energy, upregulate, and utilize PPARD. A deep understanding of its role and health and disease and how to best utilize its functions as a clinical tool is thus critical.
Species:	Human
Construct:	PPARD (His-2-441(end))
Concentration:	0.97 mg/ml
Expression System:	<i>E. coli</i>
Purity:	≥90%
Format:	Aqueous buffer solution.
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
MW:	51 kDa
Genbank Accession:	NM_006238.5
Stability:	At least 6 months at -80°C.
Storage:	-80°C
Instructions for Use:	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
Applications:	Useful for SDS-PAGE.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining

