

Recombinant PHD1 (EGLN2) protein

Catalog No: 81064, 81764

Lot No: 24717001

Expressed In: Baculovirus

Quantity: 20, 1000 µg

Concentration: 0.25 µg/µl

Source: Human

Buffer Contents: Recombinant PHD1 / EGLN2 protein is in 25 mM HEPES-NaOH pH 7.5, 300 mM NaCl, 10% glycerol, 0.04% Triton X-100 and 0.5 mM TCEP.

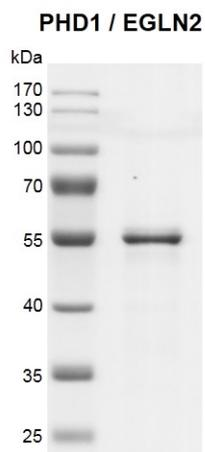
Background: Prolyl Hydroxylase Domain-Containing Protein 1 (PHD1) or Egl-9 Family Hypoxia Inducible Factor 2 (EGLN2), also known as HIF-PH1 or EIT6, is a prolyl hydroxylase. PHD3 is a cellular oxygen sensor that catalyzes the post-translational formation of 4-hydroxyproline in hypoxia-inducible factor (HIF) alpha proteins under normoxic conditions. It can hydroxylate a specific proline found in each of the oxygen-dependent degradation (ODD) domains (N-terminal, NODD, and C-terminal, CODD) of HIF1A, also HIF2A. PHD1 has a preference for the CODD site for both HIF1A and HIF2A. Hydroxylated HIFs are targeted for proteasomal degradation via the von Hippel-Lindau ubiquitination complex. Under hypoxic conditions, the hydroxylation reaction is attenuated allowing HIFs to escape degradation resulting in their translocation to the nucleus, heterodimerization with HIF1B, and increased expression of hypoxia-inducible genes. PHD1 is involved in regulating hypoxia tolerance and apoptosis in cardiac and skeletal muscle. It also regulates susceptibility to normoxic oxidative neuronal death. PHD1 links oxygen sensing to cell cycle and primary cilia formation by hydroxylating the critical centrosome component CEP192 which promotes its ubiquitination and subsequent proteasomal degradation. Besides, PHD1 can hydroxylate IKBKB, mediating NF-kappaB activation in hypoxic conditions. Its target proteins are preferentially recognized via a LXXLAP motif.

Protein Details: Recombinant human PHD1 / EGLN2 protein was expressed in a baculovirus expression system as the full length protein (accession number NP_542770.2) with an N-terminal FLAG tag. The molecular weight of the protein is 44.9 kDa.

Application Notes: This protein is suitable for use in binding assays, inhibitor screening, and selectivity profiling.

Assay Conditions: 3 µM HIF1A (HIF-1α) peptide (DLDLEALAPYIPADDDFQL) was incubated with 300 nM PHD1 protein in 30 µl reaction system containing 20 mM Tris-HCl pH 7.5, 5 mM KCl, 1.5 mM MgCl₂, 1 mM DTT, 100 µM 2-oxoglutarate, 100 µM ascorbate and 50 µM (NH₄)₂Fe(SO₄)₂·6H₂O for 2 hours at 30°C. MALDI-TOF was used for detection.

Storage and Guarantee: Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.

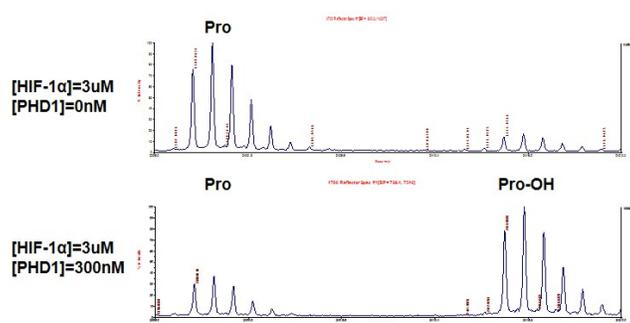


Recombinant PHD1 / EGLN2 protein gel

10% SDS-PAGE Coomassie staining

MW: 44.9 kDa

Purity: >90%



MALDI-TOF for Recombinant PHD1 / EGLN2 protein

3 μ M HIF1- α peptide was incubated with 300 nM PHD1 protein in 30 μ l reaction system for 2 hours at 30°C. The reaction product was detected by MALDI-TOF. Single 3 μ M HIF1- α peptide was used as a negative control.