

# PRODUCT INFORMATION



## PEDF (human, recombinant)

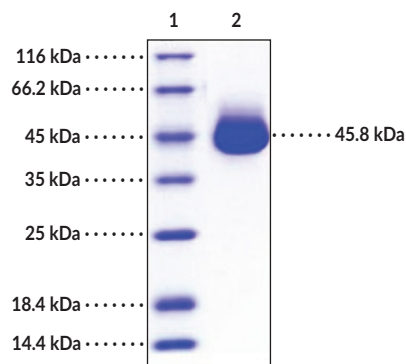
Item No. 32062

### Overview and Properties

**Synonyms:** Cell Proliferation-Inducing Gene 35 Protein, Early Population Doubling Level cDNA-1, EPC-1, Pigment Epithelium-Derived Factor, Serpin F1  
**Source:** Recombinant human C-terminal His-tagged PEDF expressed in HEK293 cells  
**Amino Acids:** 20-418  
**Uniprot No.:** P36955  
**Molecular Weight:** 45.8 kDa  
**Storage:** -80°C (as supplied)  
**Stability:** ≥1 year  
**Purity:** ≥98% estimated by SDS-PAGE  
**Supplied in:** Lyophilized from sterile PBS, pH 7.4  
**Endotoxin Testing:** <1.0 EU/μg, determined by the LAL endotoxin assay

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Image



Lane 1: MW Markers  
Lane 2: PEDF

**SDS-PAGE Analysis of PEDF.** This protein has a calculated molecular weight of 45.8 kDa.

**WARNING**  
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

**SAFETY DATA**  
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

**WARRANTY AND LIMITATION OF REMEDY**  
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 06/24/2020

**CAYMAN CHEMICAL**  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM

# PRODUCT INFORMATION



## Description

---

Pigment epithelium-derived factor (PEDF) is a secreted glycoprotein encoded by the *SERPINF1* gene.<sup>1</sup> It is a non-inhibitory member of the serine-protease inhibitor (SERPIN) superfamily that has anti-angiogenic and neurotrophic activities. PEDF is a 418-amino acid protein comprised of an N-terminal region important for its neurotrophic activity, three  $\beta$ -sheets and ten  $\alpha$ -helices that make up an  $\alpha/\beta$  core serpin domain, and a C-terminal region containing a glycosylation site.<sup>1,2</sup> It has an asymmetrical charge distribution with the basic region involved in glycosaminoglycan binding and the acidic region binding to collagen I.<sup>2</sup> PEDF is ubiquitously expressed and localized to the extracellular matrix but can also be found in the cytoplasm and nucleus of retinal pigment epithelial cells and certain cancer cell lines.<sup>1,3</sup> It contains biologically active peptides, including a 34-amino acid peptide (amino acids 44-77), which contains a laminin receptor binding site and has anti-angiogenic, bone mineralization, and apoptotic activities, and a 44-amino acid peptide (amino acids 78-121), which contains a PNPLA2 binding site and promotes neuronal differentiation and inhibits vascular permeability.<sup>1,4,5</sup> PEDF also interacts with collagen I for its anti-angiogenic activity, as well as with sulfated (heparin) and non-sulfated (hyaluronan) glycosaminoglycans, which are important for PNPLA2 receptor binding and activation of caspases, respectively.<sup>1,2</sup> Its expression is reduced in a variety of cancer cell lines and solid tumors, and the reduction is associated with metastasis and poor prognosis.<sup>2</sup> Mutations in *SERPINF1* are responsible for the hereditary bone dysplasia disorder osteogenesis imperfecta type VI, which is characterized by bone fragility, deformity, and growth deficiency.<sup>6</sup> Cayman's PEDF (human, recombinant) protein consists of 410 amino acids, has a calculated molecular weight of 45.8 kDa, and a predicted N-terminus of Gln20 after signal peptide cleavage.

## References

---

1. Kawaguchi, T., Yamagishi, S.-I., and Sata, M. Structure-function relationships of PEDF. *Curr. Mol. Med.* **10(3)**, 302-311 (2010).
2. Filleur, S., Nelius, T., de Riese, W., et al. Characterization of PEDF: A multi-functional serpin family protein. *J. Cell. Biochem.* **106(5)**, 769-775 (2009).
3. Tombran-Tink, J., Aparicio, S., Xu, X., et al. PEDF and the serpins: Phylogeny, sequence conservation, and functional domains. *J. Struct. Biol.* **151(2)**, 130-150 (2005).
4. Belinsky, G.S., Sreekumar, B., Andrejcsk, J.W., et al. Pigment epithelium-derived factor restoration increases bone mass and improves bone plasticity in a model of osteogenesis imperfecta type VI via Wnt3a blockade. *FASEB J.* **30(8)**, 2837-2848 (2016).
5. Bernard, A., Gao-Li, J., Franco, C.-A., et al. Laminin receptor involvement in the anti-angiogenic activity of pigment epithelium-derived factor. *J. Biol. Chem.* **284(16)**, 10480-10490 (2009).
6. Marini, J.C. and Blissett, A.R. New genes in bone development: What's new in osteogenesis imperfecta. *J. Clin. Endocrinol. Metab.* **98(8)**, 3095-3103 (2013).

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM