

# Progranulin (mouse), (recombinant)

Progranulin (PGRN) has been shown to be a widely expressed pluripotent growth factor which plays a role in processes such as development, wound repair and inflammation by activating signalling cascades that control cell cycle progression and cell motility. PGRN has also been linked to tumorigenesis. The function of PGRN in the central nervous system (CNS) is also of interest, because mutations in the PGRN gene have been identified in cases of frontotemporal degeneration (FTD).

**Manuals, SDS & CofA**

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	Working aliquots are stable for up to 3 months when stored at -20°C.
<b>Handling</b>	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	Proepithelin, PEPI, PC cell-derived growth factor
<b>Endotoxin Content</b>	<0.01EU/μg protein (LAL test).
<b>Formulation</b>	Lyophilized from a 0.2 μm filtered solution in PBS.
<b>Purity</b>	≥90% (SDS-PAGE)
<b>Reconstitution</b>	Reconstitute at 250 μg/ml in sterile PBS.
<b>Source</b>	Produced in mouse myeloma cell line, NS0-derived. Mouse progranulin (aa 18-589) is fused at the C-terminus with a His-tag.
<b>UniProt ID</b>	P28798

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)