

# Progranulin (human), (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).

## Ordering Information

[Order Online »](#)

ALX-201-502-0050	50µg
------------------	------

## 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>Activity</b>	Measured by its ability to enhance neurite outgrowth of E16E18 rat embryonic cortical neurons. rhProgranulin, immobilized at 1040 µg/ml on a nitrocellulose coated plate, is able to significantly enhance neurite outgrowth.
<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>MW</b>	~90-95kDa
<b>Purity</b>	≥95% (SDS-PAGE)
<b>Reconstitution</b>	Reconstitute at 200µg/ml in sterile PBS.
<b>Source</b>	Produced in mouse myeloma cell line NS0. Human progranulin (aa 18-593) is fused at the C-terminus to a 6-His tag.
<b>UniProt ID</b>	P28799

**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)