

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).

Ordering Information

[Order Online »](#)

ALX-201-503-0050	50µg
------------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

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