

Growth and Differentiation factor- 11 (human), (recombinant)

Growth differentiation factor 11 (GDF-11), is a regulator of cell growth and differentiation during muscular and neural development. GDF-11 binds the transforming growth factor-beta receptors ALK4, ALK5, and ALK7 to activate SMAD signaling.

Ordering Information

[Order Online »](#)

ENZ-PRT236-0020	20µg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -80°C. For long term storage, it is recommended to dilute to working aliquots in a 0.1% BSA solution.
Long Term Storage	-80°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	GDF-11, Bone morphogenetic protein 11, BMP-11
Biological Activity	Alkaline phosphatase activity in ATDC5 cells; ED50 \leq 100 ng/mL ($\geq 1.0 \times 10^4$ units/mg).
Endotoxin Content	≤ 1 EU/ μ g protein (LAL test)
Formulation	Lyophilized from a sterile (0.2 micron) filtered solution containing 0.1% Trifluoroacetic Acid (TFA).
MW	Dimer, $\sim 12.5/24.9$ kDa
Purity	$\geq 95\%$ (SDS-PAGE)
Reconstitution	Reconstitute in sterile 10mM HCl at 0.1 mg/mL.
Source	Produced in <i>E. coli</i> . Contains 109/218 amino acids.
UniProt ID	O95390

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)