

GDNF (human), (recombinant)

GDNF is a small protein that potently promotes the survival of many types of neurons. The most prominent feature of GDNF is its ability to support the survival of dopaminergic and motoneurons. These neuronal populations die in the course of Parkinson's disease and amyotrophic lateral sclerosis (ALS). GDNF also regulates kidney development and spermatogenesis, and it affects alcohol consumption.

Ordering Information

[Order Online »](#)

ENZ-PRT145-0010	10µg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Short Term Storage -20°C

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Glial cell-derived neurotrophic factor, Astrocyte-derived trophic factor, ATF
Formulation	Lyophilized from 10 mM sodium citrate, containing 150mM NaCl.
Purity	≥98% (SDS-PAGE gel & HPLC)
Source	Produced in <i>E.coli</i> .
UniProt ID	P39905

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)