

Neuregulin 1- β (human), (recombinant)

Neuregulin 1- β (NRG1) is one of many NRG isoforms that contains a soluble EGF-like domain. The EGF-like domain signals through receptors ErbB2, ErbB3, and ErbB4 to act as a growth factor. NRG isoforms are important for nervous system and cardiovascular development.

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.
Short Term Storage	+4°C
Long Term Storage	-80°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	NGR β 1, HRG1 β , Heregulin β 1
Endotoxin Content	≤ 0.05 EU/ μ g protein (LAL method)
Formulation	Lyophilized. Contains 0.1% trifluoroacetic acid.
MW	~ 7.6 kDa
Purity	$\geq 95\%$ (SDS-PAGE)
Reconstitution	Reconstitute in sterile water at 0.1 mg/mL.
Source	Produced in <i>E. coli</i> . A non-glycosylated protein containing 66 amino acids.
UniProt ID	Q02297-6

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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