

PRODUCT INFORMATION

Neuregulin-1 Type I-α EGF Domain (human, recombinant)

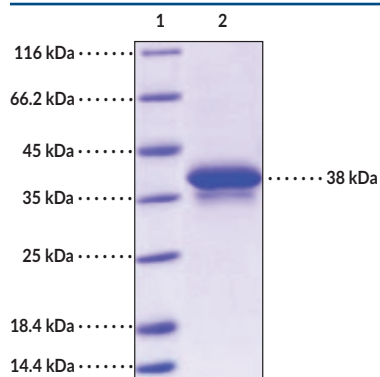
Item No. 42760

Overview and Properties

Synonyms:	Heregulin α, HRG-α, NDF-α, Neu Differentiation Factor α, NRG1-α, NRG1 Type I Isoform α
Source:	Active recombinant N-terminal human IgG1 Fc-tagged neuregulin-1 type I-α EGF domain expressed in HEK293 cells
Amino Acids:	177-241
Uniprot No.:	Q02297-1
Molecular Weight:	35.8 kDa
Storage:	-80°C (as supplied)
Stability:	≥1 year
Purity:	≥97% estimated by SDS-PAGE
Supplied in:	Lyophilized from sterile PBS, pH 7.4
Endotoxin Testing:	<1.0 EU/μg, determined by the LAL endotoxin assay
Specific Activity:	1. Measured by its ability to bind to biotinylated human HER4 in a functional ELISA. 2. Measured by its ability to bind to biotinylated rhesus monkey HER3 in a functional ELISA.

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Image



Lane 1: MW Markers
Lane 2: Neuregulin-1 Type I-α EGF Domain

SDS-PAGE Analysis of Neuregulin-1 Type I-α EGF Domain. This protein has a calculated molecular weight of 35.8 kDa. It has an apparent molecular weight of approximately 38 kDa by SDS-PAGE under reducing conditions.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 05/12/2025

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Neuregulin-1 (NRG1) type I- α is a trophic factor and member of the neuregulin-1 family of growth factors, which comprises greater than 30 isoforms grouped into types I through VI based on their N-terminal region, which is generated by alternative splicing of NRG1.¹ It is composed of a type I N-terminal domain, an immunoglobulin-like (Ig-like) domain, a glycosylated domain, an EGF-like domain, a linker region that contains a proteolytic cleavage site to release the mature protein, a transmembrane domain, and a cytoplasmic tail.^{1,2} NRG1 type I- α is expressed in brain, liver, and cardiac microvascular endothelial cells and plays a role in neural development, nerve regeneration, and gluconeogenesis.^{1,3-6} NRG1 signaling is mediated by homodimers of HER4 and the heterodimers HER2-HER3, HER2-HER4, HER3-HER4, and EGFR-HER4.¹ Low tumor levels of NRG1 type I- α are associated with decreased progression-free and overall survival in patients with breast cancer.⁷ Cayman's Neuregulin-1 Type I- α EGF Domain (human, recombinant) protein can be used for binding assays. This protein is a disulfide-linked homodimer. The reduced monomer, composed of NRG1 type I- α (amino acids 177-241) fused to human IgG1 Fc at its N-terminus, consists of 326 amino acids and has a calculated molecular weight of 35.8 kDa. The monomer migrates at approximately 38 kDa by SDS-PAGE under reducing conditions.

References

1. Mei, L. and Xiong, W.-C. Neuregulin 1 in neural development, synaptic plasticity and schizophrenia. *Nat. Rev. Neurosci.* **9**(6), 437-452 (2008).
2. Grimm, S. and Leder, P. An apoptosis-inducing isoform of neu differentiation factor (NDF) identified using a novel screen for dominant, apoptosis-inducing genes. *J. Exp. Med.* **185**(6), 1137-1142 (1997).
3. Kerber, G., Streif, R., Schwaiger, F.-W., et al. Neuregulin-1 isoforms are differentially expressed in the intact and regenerating adult rat nervous system. *J. Mol. Neurosci.* **21**(2), 149-165 (2003).
4. Arai, T., Ono, Y., Arimura, Y., et al. Type I neuregulin1 α is a novel local mediator to suppress hepatic gluconeogenesis in mice. *Sci. Rep.* **7**, 42959 (2017).
5. Cote, E.M., Miller, T.A., Lebrasseur, N.K., et al. Neuregulin-1 α and β isoform expression in cardiac microvascular endothelial cells and function in cardiac myocytes in vitro. *Exp. Cell Res.* **11**(1), 135-146 (2005).
6. De Souza, F.I., Zumiotti, A.V., and Da Silva, C.F. Neuregulins 1- α and 1- β on the regeneration the peripheral nerves. *Acta Ortop Bras.* **18**(5), 250-254 (2010).
7. Raj, E.H., Skinner, A., Mahji, U., et al. Neuregulin 1-alpha expression in locally advanced breast cancer. *Breast* **10**(1), 41-45 (2001).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM