

PRODUCT INFORMATION



Contactin-1 (human, recombinant)

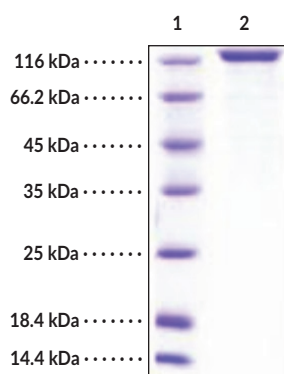
Item No. 43685

Overview and Properties

Synonyms: CNTN1, GP135, Neural Cell Surface Protein F3
Source: Recombinant human C-terminal His-tagged contactin-1 expressed in HEK293 cells
Amino Acids: 21-993
Uniprot No.: Q12860
Molecular Weight: 110 kDa
Storage: -80°C (as supplied)
Stability: ≥1 year
Purity: ≥95% 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
Protein Concentration: *batch specific* mg/ml

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

Image



Lane 1: MW Markers
Lane 2: Contactin-1

SDS-PAGE Analysis of Contactin-1. This protein has a calculated molecular weight of 110 kDa. It has an apparent molecular weight of approximately 125 kDa due to glycosylation.

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, 06/11/2025

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

Contactin-1 is a neural cell adhesion molecule and member of the immunoglobulin (Ig) superfamily.¹ It is composed of six Ig-like repeats, four fibronectin type III-like domains, and a hydrophobic C-terminal sequence that contains a glycosylphosphatidylinositol (GPI) anchor site.^{1,2} Contactin-1 is primarily expressed in the brain and to a lesser extent in lung, pancreas, skeletal muscle, and kidney.² It is involved in many neuronal developmental processes, including axonal and neurite growth, glial cell differentiation, myelination, and synaptogenesis.^{1,3} Contactin-1 has several interaction partners, such as Notch, protein tyrosine phosphatase α (PTP α), and tenascin-R, and binds to them in either *cis* or *trans*.³ Increased expression of *CNTN1*, the gene encoding contactin-1, is associated with decreased overall survival in patients with bladder urothelial carcinoma or stomach adenocarcinoma.¹ Mutations in *CNTN1* are associated with familial lethal congenital myopathy.⁴ Cerebrospinal fluid (CSF) levels of contactin-1 are decreased in patients with Parkinson's disease.⁵ Cayman's Contactin-1 (human, recombinant) protein consists of 984 amino acids, has a calculated molecular weight of 110 kDa, and a predicted N-terminus of Glu21 after signal peptide cleavage. By SDS-PAGE, under reducing conditions, the apparent molecular mass of the protein is 125 kDa due to glycosylation.

References

1. Gu, Y., Li, T., Kapoor, A., *et al.* Contactin 1: An important and emerging oncogenic protein promoting cancer progression and metastasis. *Genes (Basel)* **11(8)**, 874 (2020).
2. Reid, R.A., Bronson, D.D., Young, K.M., *et al.* Identification and characterization of the human cell adhesion molecule contactin. *Brain Res. Mol. Brain Res.* **21(1-2)**, 1-8 (1994).
3. Shimoda, Y. and Watanabe, K. Contactins: Emerging key roles in the development and function of the nervous system. *Cell Adh. Migr.* **3(1)**, 64-70 (2009).
4. Compton, A.G., Albrecht, D.E., Seto, J.T., *et al.* Mutations in contactin-1, a neural adhesion and neuromuscular junction protein, cause a familial form of lethal congenital myopathy. *Am. J. Hum. Genet.* **83(6)**, 714-724 (2008).
5. Chatterjee, M., van Steenoven, I., Huisman, E., *et al.* Contactin-1 is reduced in cerebrospinal fluid of Parkinson's disease patients and is present within Lewy bodies. *Comparative Study* **10(8)**, 1177 (2020).

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